•TECAN•

Instrucciones de uso de



magellan Tracker:	7.5
magellan:	7.5
Revisiones del software:	
Nº de versión del documento:	1.5
	2023-07
Nº de pieza del documento:	30143535

CE



30143535 05



ADVERTENCIA ANTES DE UTILIZAR EL INSTRUMENTO, LEA CON CUIDADO EN ESTE MANUAL Y APLIQUE LAS INSTRUCCIONES DADAS.

Aviso

Hemos realizado el máximo esfuerzo para evitar errores en el texto y en los diagramas; no obstante, Tecan Austria GmbH no asume ninguna responsabilidad por posibles errores en la presente publicación.

Tecan Austria GmbH tiene el compromiso de mejorar nuestros productos para aprovechar las nuevas técnicas y componentes. Por eso nos reservamos el derecho de cambiar las especificaciones en cualquier momento, cumpliendo las correspondientes verificaciones, validaciones y homologaciones.

Agradecemos recibir comentarios sobre la presente publicación.



Fabricante

Tecan Austria GmbH Untersbergstr. 1A A-5082 Grödig, Austria T: +43 6246 89330 www.tecan.com E-mail: office.austria@tecan.com

Información de copyright

El contenido de este manual es propiedad de Tecan Austria GmbH y no se puede copiar, reproducir ni transferir a terceros sin nuestro consentimiento previo por escrito.

Copyright © Tecan Austria GmbH Reservados todos los derechos. Impreso en Austria.

Declaración de certificado UE

Se suministra a petición.

Uso especificado de magellan

Consulte 1.2 Uso / fin especificado de magellan.

Acerca de este manual

magellan es un paquete universal de reducción de datos que se utiliza para analizar los datos generados a partir de ensayos con microplacas. Está diseñado para uso profesional.

Este manual explica cómo:

- Instalar el software
- Utilizar el software

Nota acerca de las capturas de pantalla

Es posible que el número de versión que muestran algunas capturas de pantalla no siempre coincida con el de la versión más actualizada. Las capturas de pantalla solamente estarán actualizadas si ha cambiado el contenido correspondiente de la aplicación.



Historial de cambios de las Instrucciones de utilización

Versión	Cambios
V 1.2	Nueva DoC (Declaration of Conformity)

Advertencias, Precauciones y Avisos

En esta publicación se usan los tres tipos de aviso siguientes para hacer resaltar información importante o alertar al usuario sobre situaciones potencialmente peligrosas:





Símbolos

	Fabricante
\sim	Fecha de fabricación
CE	Marcado CE de conformidad
ĺ	Lea las Instrucciones de uso antes de manejar este instrumento
IVD	Producto sanitario para diagnóstico in vitro
UK CA	United Kingdom Conformity Assessed El marcado "United Kingdom Conformity Assessed" muestra que el producto etiquetado cumple la reglamentación aplicable en Gran Bretaña.
UDI	Unique Device Identification El símbolo UDI (Identificación única de dispositivos) indica al portador de datos en la etiqueta.
REF	Número de pedido
SN	Número de serie
X	Símbolo de WEEE
	Símbolo de RoHS, China



Contenido

1.	Introducción9			
	1.1	Campo) de aplicación	9
	1.2	Uso / fi	in especificado de magellan	10
	1.3	Perfiles	s de usuarios	10
		1.3.1	Usuario profesional - Nivel Administrador	10
		1.3.2	Usuario final o usuario rutinario	10
	1.4	Especit	ficaciones	
		1.4.1	Requisitos del sistema	11
		1.4.2	Compatibilidad con lectores	11
	1.5	Proced	limiento de instalación del software	
		1.5.1	Programa de instalación automática del software	12
		1.5.2	Comprobación de la instalación – IQ	
		1.5.3	Validación de operatividad (OQ, Operational Qualification)	
		1.5.4	Recuperación del sistema	
		1.5.5	Eliminación automática del software	
		1.5.6	Información sobre seguridad cibernética	
2.	Emp	oezar a tr	rabajar con magellan	17
	2.1	Interfaz	z del usuario – Lista de asistentes	17
	2.2	Compo	onentes y términos – Lógica básica de magellan	
		2.2.1	Tipos de archivo usados con magellan	
		2.2.2	Procesamiento de carpetas	
		2.2.3	Elementos estándar	21
		2.2.4	El botón de Ayuda	
		2.2.5	El cuadro de diálogo de bienvenida	
		2.2.6	Lista de accesos directos	
	2.3	Iniciar	magellan	
		2.3.1	Iniciar la versión estándar	
		2.3.2	Iniciar la versión Tracker	
	2.4	Conect	tar un instrumento	
		2.4.1	Conectarse a instrumentos SUNRISE	
		2.4.2	Conectarse a instrumentos INFINITE F50	
		2.4.3	Registrar magellan	
	_	2.4.4	Asistente para registrar	
3.	Con	trol y co	nfiguración del instrumento	31
	3.1	Opcion	nes del instrumento	
		3.1.1	Movimientos	
		3.1.2	Control de temperatura	
	3.2	Opcion	nes de Configuración y Servicios	
		3.2.1	Cambiar instrumento	
		3.2.2	Definir placas de filtro	
	3.3	Archive	os de registro	
4.	Asis	stente Cr	rear o editar método	33
	4.1	Introdu	icción	
	4.2	Definir	los parámetros de medición	
		4.2.1	Tipos de mediciones - SUNRISE	
		4.2.2	Parámetros de medición - SUNRISE	
		4.2.3	Parámetros de medición - INFINITE F50	
	4.3	Definir	evaluación	39
		4.3.1	Ventana de visión global de Crear o editar método	
		4.3.2	Diseño de método: Cómo definir un diseño de placas	
		4.3.3	Diseño de método: Valores Conc., Dil. y Ref	
		4.3.4	Cálculo previo: Reducción de datos de espectros	50



		4.3.5	Datos transformados: Añadir nueva transformación	53
		4.3.6	Cinética: Reducción de datos cinéticos	56
		4.3.7	Transformaciones cinéticas: Añadir nuevas transformaciones cinéticas	60
		4.3.8	Concentraciones: Curva estándar	60
		4.3.9	Transformaciones de concentraciones: Añadir nueva transformación de concentración.	66
		4.3.10	Evaluar datos: Definición del corte	66
		4.3.11	Evaluar datos: Validación QC	68
		4.3.12	Procesamiento de datos: Exportación de datos	69
		4.3.13	Procesamiento de datos: Informe impreso	74
		4.3.14	Procesamiento de datos: Procesamiento automatizado de datos	79
		4.3.15	Otras opciones: Indicaciones al usuario	84
		4.3.16	Otras opciones: Formato de números	84
		4.3.17	Otras opciones: Notas de método	85
	4.4	Guardar	[•] métodos	85
		4.4.1	Protección de métodos mediante contraseña	86
	4.5	Métodos	s multiplaca	87
5.	Asis	tente Cre	ear o editar lista de IDs de muestras	89
	5.1	Introduc	ción	89
	5.2	Crear o	editar lista de IDs de muestras	89
		5.2.1	Crear nueva lista de IDs de muestras	89
		5.2.2	Importar o modificar lista de IDs de muestras	91
		5.2.3	Importar lista de IDs de muestras	96
		5.2.4	Guardar listas de IDs de muestras	104
6.	Asis	tente Inic	ciar medición	.105
	6.1	Introduc	ción	105
	6.2	Obtener	datos sin procesar	106
		6.2.1	Obtener datos sin procesar con el instrumento SUNRISE	106
		6.2.2	Obtener datos sin procesar con el instrumento INFINITE F50	106
	6.3	Usar mé	etodo predefinido	107
	6.4	Activar	favorito	107
	6.5	Iniciar m	nedición con un método predefinido o favorito	108
-	6.6	Estado o	de la medición	
1.	ASIS	tente Eva	aluar resultados	.115
	7.1	Introduc	ccion	115
	7.2	Seleccio	onar un archivo	115
	7.0	7.2.1 Mantana	Criterios de selección de archivos	116
	7.3	Ventana	l vista de conjunto de la zona de trabajo	117
	7.4		Evaluar resultados	
		7.4.1	Menus	120
		7.4.2	Menú de Barra de herramientas: Alchivo	120
		7.4.3	Menú de Barra de herramientas: Eulcion	124
		7.4.4	Venteno de diseño de places	125
		7.4.5		125
		7.4.0	Calacteres especiales	120
		1.4.1 710	Barra de control: Datos reducidos	120 196
		7.4.0	Barra de control: Datos transformados	120
		7 <u>1</u> 10	Barra de control: Parámetros cinéticos	120
		7 4 11	Barra de control: Concentraciones	120
		7 4 12	Barra de control: Resultados cualitativos	1.21
		7412	Barra de control: IDs de muestras	1.31
		7.4.14	Barra de control: Diseño de método	
		7.4.15	Barra de control: Validación QC	132
		7.4.16	Barra de control: Otras opciones	132



		7.4.17	Cuadro de diálogo Escala de colores	132
		7.4.18	Menú contextual de un pocillo	132
	7.5	Pestaña	Editar método	137
	7.6	Guardar	los Resultados evaluados	138
8.	Asist	tente Ane	exar firma	139
	8.1	Introduc	ción	139
	8.2	Firmar u	n archivo	139
9.	Icono	o Otras o	pciones	141
	9.1	Control	del instrumento	
	9.2	Procesa	miento de archivos	141
		9.2.1	Guardar en archivador	141
		9.2.2	Importar datos sin procesar	144
		9.2.3	Convertir a	145
		9.2.4	Convertir de	146
		9.2.5	Guardar archivos de registro	147
	9.3	Opcione	S	
		9.3.1	Directorios predeterminados de datos	
		9.3.2	Opciones de Copiar/Exportar	150
		9.3.3	Pestaña Vista de placa	152
		9.3.4	Otras opciones	153
	9.4	Adminis	tración de usuarios (magellan Tracker)	155
		9.4.1	Añadir/modificar usuario (magellan Tracker)	156
		9.4.2	Agregar/modificar grupo	
		9.4.3	Registro de modificaciones de administración de usuarios	
		9.4.4	Opciones de Administración de usuarios	
		9.4.5	Resumen de Administración de usuarios	
	9.5	Adminis	tración de usuarios (magellan Standard)	
		9.5.1	Anadir/modificar usuario (magelian Standard)	
		9.5.2	Inicio de sesion	
	0.0	9.5.3 Dereche	Campiar usuario	
	9.0	Derecho	s dei usuario	
10	9.7 Euno	Acerca d	de magellan	
10.	Func 10.1			
	10.1	Auminis	Pagiatra da madificacianas	171
	10.2	Drocces	Registro de modificaciones	
	10.2	10 2 1	Guardar un archivo	
		10.2.1	Cambiar un archivo de zona de trabajo o mátodo	
		10.2.2	Abrir un archivo	
		10.2.3	Abrir un archivo creado en otro ordenador – Añadir HIIDs	
	10 3	Registro	de modificaciones del sistema	173
11.	Cálci			
• • •	11.1	Fvaluar	resultados – Procedimiento de cálculo	175
		11.1.1	Estadísticas	
	11.2	Reducci	ón de datos de espectros	
		11.2.1	Descripción matemática	
	11.3	Escribir	una fórmula	
	-	11.3.1	Introducción	
		11.3.2	Variables en fórmulas	
		11.3.3	Funciones de fórmulas	
		11.3.4	Funciones básicas	
		11.3.5	Funciones estadísticas	
		11.3.6	Funciones de eliminación	
		11.3.7	Otras funciones	187



		11.3.8	Funciones de espectros	188
		11.3.9	Ejemplos	189
	11.4	Tipos de	análisis de la curva estándar	190
		11.4.1	Definiciones	190
		11.4.2	Parámetros de tipo de análisis	190
		11.4.3	Mensajes de error	191
		11.4.4	Punto a punto	191
		11.4.5	Regresión lineal	191
		11.4.6	Regresión no lineal	192
		11.4.7	Polinomio	193
		11.4.8	Spline cúbica	194
		11.4.9	Akima	194
		11.4.10	LogitLog	195
		11.4.11	Cuatro parámetros	197
		11.4.12	Cuatro parámetros – Marquardt	197
		11.4.13	Cinco parámetros – Marquardt	198
		11.4.14	Ponderación para ajuste de cuatro / cinco parámetros – ajuste polinomial / Marquardt	199
	11.5	Cálculo d	de series de dilución	200
		11.5.1	Detección de series de dilución	200
		11.5.2	Cálculo de parámetros de curvas	200
		11.5.3	Cálculo de valores de IC	200
12.	Ejem	plo de ap	olicación	201
	12.1	Ejemplo	paso a paso: ELISA cuantitativo	201
		12.1.1	Descripción del kit de prueba	201
		12.1.2	Crear un método	202
		12.1.3	Ejecutar el método	216
		12.1.4	Evaluar el resultado	217
		12.1.5	Resumen de la definición de ELISA cuantitativo en magellan	219
13.	Glos	ario		221
Índic	e			225
Marc	as reg	jistradas.		228
Aten	ción a	I Cliente	de Tecan	229



1. Introducción

1.1 Campo de aplicación

El software **magellan** es un programa de control de lector y análisis de datos para analizar datos generados por mediciones realizadas con un lector de absorbencia de microplacas monomodo de Tecan (por ejemplo, SUNRISE e INFINITE F50) para diagnósticos in vitro.



Aviso Es importante tener en cuenta que la correcta instalación de este instrumento y el software magellan no garantiza en sí misma la conformidad con todas las leyes y requisitos legales. También es preciso establecer las medidas correspondientes respecto a los procesos y procedimientos operativos estándar, que incluyen la validación y el control de calidad.

magellan está diseñado para usarse con las versiones de sistema operativo y Microsoft Office especificadas en el capítulo 1.4.1 Requisitos del sistema. Si se instalan programas adicionales, no se puede garantizar la funcionalidad según las Regulaciones IVD para Europa y FDA.

Aviso

Para evitar el uso indebido de derechos de usuario y la falsificación de datos, es recomendable que el Administrador de usuarios no tenga derechos en Magellan. A ser posible, el Administrador de usuarios deberá formar parte del departamento de TI.

Aviso

El administrador del sistema de los clientes es responsable de los cambios realizados en el sistema operativo de los equipos. El cliente tiene que asegurarse de establecer los derechos de usuario apropiados para evitar la eliminación o la manipulación de datos fuera del software magellan.

Aviso magellan está diseñado para el uso en un solo ordenador, no integrado en una red.

i

Aviso

Cualquier incidente grave sucedido en relación con el dispositivo y el software se notificará al fabricante y a la autoridad competente del Estado miembro donde se encuentre establecido el usuario o el paciente.

1.2 Uso / fin especificado de magellan

El software **magellan** es un programa de control de lector y análisis de datos para analizar datos generados por mediciones realizadas con un lector de absorbencia de microplacas monomodo de Tecan (por ejemplo, SUNRISE e INFINITE F50) para diagnósticos in vitro.

El software **magellan** está destinado principalmente a mediciones de absorbencia y la evaluación de ensayos IVD cuantitativos, semicuantitativos y cualitativos. El software **magellan** es programable por el usuario, que debe configurarlo de acuerdo con las instrucciones de cada ensayo.

Este producto está destinado al uso profesional en laboratorio por personal capacitado, no es para uso doméstico ni por parte de profanos.

magellan está disponible en dos versiones:

- magellan
- magellan Tracker

magellan Tracker ofrece toda la funcionalidad necesaria para cumplir el Reglamento de la FDA 21 CFR parte 11, además de la funcionalidad de **magellan**.



Aviso magellan está disponible en dos versiones: magellan y magellan Tracker. Para facilitar la referencia en este manual de instrucciones, la versión de

magellan se marca como magellan Standard.

1.3 Perfiles de usuarios

1.3.1 Usuario profesional - Nivel Administrador

El administrador es una persona que tiene la formación técnica apropiada y las cualificaciones y la experiencia correspondientes. Si el equipo se utiliza como está previsto, esa persona podrá reconocer y evitar el peligro.

El Administrador deberá tener amplios conocimientos y podrá instruir al usuario final y al usuario rutinario sobre los protocolos de análisis en relación con un producto de Tecan dentro de los límites del uso previsto.

Se requieren destrezas informáticas y el idioma inglés.

1.3.2 Usuario final o usuario rutinario

El usuario final (o usuario rutinario) es una persona que tiene la formación técnica apropiada y las cualificaciones y la experiencia correspondientes. Si el equipo se utiliza como está previsto, esa persona podrá reconocer y evitar el peligro.

Se requieren destrezas informáticas y buenos conocimientos del idioma local respectivo o inglés.



Aviso Consulte en Atención al cliente los detalles sobre formación (fechas, duración y frecuencia). Puede encontrar las direcciones y números de teléfono en Internet: http://www.tecan.com/customersupport

1.4 Especificaciones

1.4.1 Requisitos del sistema

Para utilizar magellan se deben cumplir los siguientes requisitos:

	Mínimo	Recomendado
PC	PC compatible con Windows y procesador compatible con Pentium a 1 GHz	2 GHz (Doble núcleo)
Sistema operativo	Windows 10 (32 bits)	
	Windows 10 (64 bits): Ediciones: Pro	
Memoria	Windows 10 (32 bits): 1 GB de RAM	2 GB de RAM
	Windows 10 (64 bits): 2 GB de RAM	4 GB de RAM
Espacio libre en el disco duro	3 GB	5 GB
Monitor	Gráficos SuperVGA	
Resolución	1024 x 600 y posterior	1920 x 1080
Profundidad de color	256	
Ratón	Microsoft Mouse o dispositivo señalador compatible mejorado	
Comunicaciones	1 x USB 2.0	2 x USB 2.0 1 x RS232 (Serie)
Dispositivos	Windows 10: Dispositivo de gráficos DirectX 9 con controlador WDDM 1.0 o posterior	
.NET	Microsoft .NET Framework 3.5: En Windows 10, si todavía no está instalada, se pedirá al usuario que instale la versión de .NET Framework requerida (3.5).	
Windows Installer	3.1 Si esta versión no está presente, la instalará el programa de instalación o actualización.	
Microsoft Excel	2007, 2010, 2013, 2016 (32 bits), 2019 (32 bits), Microsoft Excel 365 (32 bits) ¡Solo son compatibles las ediciones de 32 bits! ¡Las ediciones Starter NO son compatibles!	2010 (32 bits) 2019 (32 bits)

1.4.2 Compatibilidad con lectores

Se pueden usar los lectores Tecan siguientes al utilizar magellan:

Tipos de instrumentos	Modos de medición
SUNRISE	Absorbencia
INFINITE F50	Absorbencia

1.5 Procedimiento de instalación del software

1.5.1 Programa de instalación automática del software

magellan se instala mediante un programa de instalación automática de software. Todos los componentes necesarios se instalan automáticamente.

Ejecute Magellan_Setup.exe para iniciar el procedimiento de instalación:

- En la primera ventana del Asistente InstallShield se muestran los componentes que es necesario instalar antes de instalar Magellan. Haga clic en Install [Instalar] para continuar.
 Windows 10: Si no está presente la versión de .NET Framework (3.5) requerida, se pide al usuario que inicie su instalación. Seleccione Install this feature [Instalar esta característica] (se requiere una conexión a Internet). Si el usuario omite esa instalación, la instalación de Magellan no será correcta.
- Se muestra una **Ventana de bienvenida** que contiene información acerca del software y el proceso de instalación. Haga clic en **Next** [Siguiente] para continuar.
- La siguiente ventana muestra el License Agreement [Contrato de licencia]. Léalo y haga clic en l agree [Acepto] para aceptarlo y continuar.
- A continuación, se muestra la ventana Customer Information [Información del cliente]. Si usted ya ha adquirido el software, introduzca ahora su número de serie y haga clic en Next. Si todavía no dispone de un número de serie, haga clic en Next igualmente para continuar con la instalación. Podrá registrarse más adelante (consulte el capítulo 2.6 Licencias de magellan).
- A continuación, se muestra la ventana Destination Folder [Carpeta de destino]. Se mostrará la ruta de instalación predeterminada. Haga clic en Browse [Examinar] para cambiar la ruta de destino predeterminada, si lo desea. Haga clic en Next para continuar.
- A continuación, se muestra la ventana Language Selection [Selección de idioma]. Seleccione su idioma preferido.
- En la ventana **Use For Regulated Environments** [Uso para entornos regulados], seleccione magellan **Standard** o magellan **Tracker**, la versión que haya solicitado, para instalar el software. Haga clic en **Next** para continuar.
- El programa de instalación ahora está listo para la instalación. Haga clic en **Next** para comenzar la instalación.
- Haga clic en **Finish** [Finalizar] para terminar la instalación y cerrar el programa de instalación.

Para iniciar el software a través del menú **Inicio** de Windows, seleccione **magellan**.

El programa de instalación detecta automáticamente una instalación previa de magellan. Antes de instalar una versión nueva de Magellan es preciso desinstalar la versión antigua. La desinstalación se puede realizar automáticamente utilizando el Asistente de instalación (ver también el capítulo 1.5.5 Eliminación automática del software).



Aviso

Es de suma importancia que la persona que instale el software tenga derechos de administrador en el ordenador.





Aviso Si desea instalar una versión anterior de magellan, previamente debe desinstalar por completo la versión actual.

1.5.2 Comprobación de la instalación – IQ

Es recomendable utilizar el programa automático de validación de la instalación para comprobar que la instalación de **magellan** ha sido correcta.

Inicie Magellan IQ.exe desde el menú Inicio de Windows: Inicio > Programas > Tecan > Magellan IQ o directamente desde el sistema de archivos utilizando la ruta de instalación predeterminada que se especificó en el programa de instalación (normalmente, C:\Program Files\Tecan\Magellan).

El programa de validación de la instalación detecta automáticamente la versión instalada de **magellan**. Haga clic en **Verificar** para iniciar la validación de la instalación.

La herramienta IQ muestra el estado de todos los componentes instalados. Hay tres estados posibles: **Correcto, Advertencia y Fallo (=Error)**. En caso de que se notifiquen errores, póngase en contacto con su departamento de soporte local para solicitar ayuda.

Después de que finalice la comprobación, se puede generar un informe que contiene la información que muestra la herramienta IQ. Además de la información de la herramienta IQ, el informe contiene un campo de firma, de modo que el informe se puede guardar e imprimir para auditoría. Haga clic en el botón **Informar** en: **Archivo > Informar** para generar un informe. El informe se puede guardar como archivo PDF y en otros formatos.

Haga clic en **Cancelar** o en **Salir** para cerrar el programa de validación de la instalación.



Aviso La validación de la instalación debe repetirse cada vez que se instale o se actualice magellan a una nueva versión, o bien se actualice o modifique el sistema subyacente.

1.5.3 Validación de operatividad (OQ, Operational Qualification)

Además de comprobar la validación de la instalación, Tecan también recomienda comprobar las capacidades de cálculo de **magellan**. Para este fin, el medio de instalación de **magellan** contiene una zona de trabajo y un archivo de informe.

Abra el archivo de zona de trabajo OQ para la versión de **magellan** instalada e imprima el informe. El informe impreso debe ser idéntico al correspondiente informe de OQ en formato PDF (por supuesto, exceptuando la información de fecha y hora).

Si esos dos informes difieren, póngase en contacto con su distribuidor local para solicitar ayuda.

1.5.4 Recuperación del sistema

Para poder reparar una instalación de **magellan** dañada después de un fallo de hardware, es importante efectuar con antelación una copia de seguridad de los datos.



Recuperación del sistema con software de copia de seguridad de datos

La copia de seguridad de los datos normalmente se lleva a cabo con un software de copia de seguridad de datos. Dicho software permite guardar el sistema completo, incluyendo todos los datos, de modo que se puedan restaurar siempre que sea necesario.

Puede encontrar una descripción detallada e instrucciones de usuario precisas en la ayuda de Windows (para paquetes de software de copia de seguridad de Windows) o en la documentación suministrada.

La licencia de **magellan** está asociada al hardware del ordenador, lo cual significa que podría invalidarse después de cambiar el hardware. Por lo tanto, después de restaurar el sistema es preciso comprobar que la licencia sigue siendo válida. Puede hacerlo en el cuadro de diálogo **Acerca de magellan**. Para ello, inicie el asistente para registrar y compruebe que el número de serie y el número de licencia registrados siguen siendo válidos. En caso de recibir un mensaje de error, póngase en contacto con el servicio de atención al cliente de Tecan para obtener un número de licencia nuevo.

Recuperación manual del sistema

En aquellos casos en los que no es posible una restauración completa del sistema con el software de copia de seguridad (p. ej., porque los cambios de hardware son demasiado numerosos o porque se ha instalado una versión distinta del sistema operativo Windows en un nuevo ordenador), es preciso reinstalar **magellan** y recuperar varios archivos desde una copia de seguridad.

En ese caso, el proceso de restauración del sistema consta de estos pasos:

- 1. Instalar magellan
- 2. Iniciar magellan y definir el Administrador de magellan
- 3. Volver a introducir el usuario y los derechos de usuario.
- 4. Copiar del archivo de copia de seguridad el archivo magellan.ini.
- 5. Copiar del archivo de copia de seguridad los métodos de magellan.
- 6. Copiar del archivo de copia de seguridad las listas de IDs de muestras.
- 7. Copiar del archivo de copia de seguridad las zonas de trabajo existentes.
- 8. Copiar del archivo de copia de seguridad los datos exportados.

Después de implementar los cambios de hardware y completar la introducción de las licencias de magellan Tracker, al abrir archivos de métodos y de zonas de trabajo generados bajo la licencia antigua se notificará al usuario que los archivos proceden de un ordenador distinto. Para solucionar este problema, siga la descripción detallada del capítulo 10.2.4 Abrir un archivo creado en otro ordenador – Añadir HUIDs.

Seguridad de los datos

Para asegurarse de que pueda completar rápidamente los pasos del proceso de restauración del sistema enumerados anteriormente, deberá realizar a tiempo una copia de seguridad. La copia de seguridad deberá incluir métodos, zonas de trabajo, listas de IDs de muestras, datos exportados, archivos de registro de modificaciones, archivos de registro de modificaciones del sistema, el archivo UserManagement.xml, archivos del directorio Pdfx y el archivo magellan.ini de **magellan**. El contenido de la copia de seguridad siempre debería estar actualizado. Tecan recomienda utilizar la opción **Autoarchivado** para hacer copias de seguridad de métodos, zonas de trabajo y listas de IDs de muestras. Puede encontrar detalles acerca del Autoarchivado en el capítulo 9.2.2 Importar datos sin procesar.



Magellan.ini	C:\Users\Public\Documents\Tecan\Magellan
Métodos	Ruta estándar: C:\Users\Public\Documents\Tecan\Magellan\mth o como se defina en el cuadro de diálogo Onciones
Zonas de trabajo	C:\Users\Public\Documents\Tecan\Magellan\wsp o como se defina en el cuadro de diálogo Opciones
Listas de IDs de muestras	C:\Users\Public\Documents\Tecan\Magellan\smp o como se defina en el cuadro de diálogo Opciones
Datos exportados	C:\Users\Public\Documents\Tecan\Magellan\asc o como se defina en el cuadro de diálogo Opciones
Archivos del registro de modificaciones	Ruta de acceso tal como se muestra en el cuadro de diálogo Registro de modificaciones (en Administración de usuarios)
Registro de modificaciones del sistema	Ruta de acceso según se especifica en 10.3 Registro de modificaciones del sistema
Administración de usuarios	C:\ProgramData\Tecan\Tecan User Management\v1.1\UserManagement.xml
Archivos del directorio Pdfx	C:\Users\Public\Documents\Tecan\Pdfx

Directorios de magellan – Windows 10

1.5.5 Eliminación automática del software

El software **magellan** puede eliminarse mediante la rutina de desinstalación estándar de Windows:

- Cierre magellan
- Seleccione Agregar o Quitar Programas desde el Panel de Control en el menú Inicio de Windows.
- Seleccione el icono de magellan y haga clic en Quitar.

Durante el proceso de eliminación se le preguntará si desea eliminar los componentes compartidos. Si selecciona **S**í, **magellan** se desinstala completamente y no funcionarán otros programas de Tecan. Si se dejan instalados esos componentes, el sistema no se verá afectado. En caso de duda, se recomienda no eliminar los componentes compartidos. Después de elegir **S**í o **No**, se desinstala el software **magellan**.



Aviso Al eliminar los componentes compartidos, también se eliminan los datos de administración de usuarios. Algunos programas de Tecan, como por ejemplo EVOware, no funcionarán más y tienen que volverse a instalar.

1.5.6 Información sobre seguridad cibernética

PRECAUCIÓN

ES RESPONSABILIDAD DEL USUARIO TOMAR LAS PRECAUCIONES NECESARIAS PARA PROTEGER EL SISTEMA FRENTE A AMENAZAS DE SEGURIDAD CIBERNÉTICA.

AL INSTALAR SOFTWARE DE PROTECCIÓN CONTRA VIRUS O ACTUALIZACIONES RELACIONADAS CON LA SEGURIDAD DEL SISTEMA OPERATIVO, SIGA LAS RECOMENDACIONES DE SU DEPARTAMENTO LOCAL DE TI. HAGA COPIA DE SEGURIDAD DE LOS ARCHIVOS DE MAGELLAN ANTES DE COMENZAR LAS ACTUALIZACIONES. DESPUÉS DE MODIFICAR EL SISTEMA, EJECUTE INSTALLATION QUALIFICATION (IQ), OPERATION QUALIFICATION (OQ) Y PERFORMANCE QUALIFICATION (PQ) DE MAGELLAN.

Haga copia de seguridad de sus datos

Es preciso insistir en la importancia vital de realizar copias de seguridad de los archivos para poder recuperarlos en caso de un ataque de ransomware. Asegúrese de crear copias de seguridad indestructibles de sus sistemas y datos críticos. Hay muchas herramientas y servicios disponibles para la copia de seguridad y la restauración de versiones anteriores los archivos, así como la recuperación ante desastres. Asegúrese también de comprobar regularmente que las copias de seguridad se están realizando.

Utilice la funcionalidad de archivado de magellan o del producto que elija para realizar copias de seguridad de los archivos de datos de magellan. Para obtener una descripción más detallada, consulte el capítulo 9.2.1 Guardar en archivador.

Administración y control del acceso con privilegios a datos

Para minimizar el riesgo de abusos y credenciales transgredidas, implemente contraseñas de administrador locales únicas en todos los sistemas, separe y proteja las cuentas con privilegios y reduzca los permisos generales de los repositorios de archivos.

Al utilizar magellan, trabaje iniciando sesión con usuarios estándar de Windows (no administradores).

Use la funcionalidad de magellan Tracker para controlar el acceso y asegurar aún más la aplicación.

Validación de operatividad

Además de verificar la validación de la instalación, Tecan también recomienda que se comprueben las capacidades de cálculo de magellan. Para obtener una descripción más detallada, consulte el capítulo 1.5.3 Validación de operatividad (OQ, Operational Qualification).



2. Empezar a trabajar con magellan

La interfaz de usuario principal en **magellan** es el asistente.

Los asistentes estándar de **magellan** representan módulos de flujo de trabajo con instrucciones paso a paso para ejecutar procedimientos complejos.

A veces los menús están disponibles en la barra de cabecera. El **Menú** ofrece una forma tradicional de usar el software: el elemento del menú correspondiente se selecciona desde los menús principales. Todas las acciones subsiguientes se iniciarán de inmediato, o bien se mostrará un cuadro de diálogo para poder seleccionar opciones adicionales.

2.1 Interfaz del usuario – Lista de asistentes

La interfaz de usuario principal en magellan es el asistente.

Después de iniciar magellan, aparece la Lista de asistentes:

\frown	¿Qué desea hacer?
Iniciar medición	 El asistente Iniciar medición le ayuda a realizar una medición. Puede utilizar un método u obtener datos
Evaluar resultados	O sin procesar
Adjuntar firma	0
Crear o editar lista de IDs de muestras	0
Crear o editar método	0
/	
Salir Magellan	

Cada asistente se puede iniciar haciendo doble clic o seleccionándolo y luego

pulsando el botón Siguiente 🕑.

Asistente Iniciar medición

El asistente Iniciar medición incluye las siguientes opciones:

- **Obtener datos sin procesar** se utiliza para establecer los parámetros de medición necesarios y generar datos en bruto de forma rápida y sencilla.
- Usar método predefinido se utiliza para realizar mediciones basadas en métodos definidos previamente.
- Activar favorito se utiliza para seleccionar uno de los métodos más frecuentes en la lista de iconos enumerados.

Para obtener una descripción más detallada, consulte el capítulo 6 Asistente Iniciar medición.

Cuando finaliza la medición, se crea un archivo de zona de trabajo.



Asistente Evaluar resultados

El **asistente Evaluar resultados** se utiliza para ver los datos en bruto y evaluar los resultados. Se pueden ver los parámetros de evaluación y pueden volverse a evaluar.

Para obtener una descripción más detallada, consulte el capítulo 7 Asistente Evaluar resultados.

Asistente Anexar firma

El **asistente Anexar firma** se utiliza para firmar archivos de método y zona de trabajo. Esta función solo está disponible con **magellan** Tracker.

Para obtener una descripción más detallada, consulte el capítulo 8 Asistente Anexar firma.

Asistente Crear o editar lista de IDs de muestras

El **asistente Crear o editar lista de IDs de muestras** se utiliza para crear listas de IDs de muestras nuevas o para editar las listas existentes.

Para obtener una descripción más detallada, consulte el capítulo 5 Asistente Crear o editar lista de IDs de muestras.

Asistente Crear o editar método

El asistente Crear o editar método se utiliza para definir o editar métodos.

Para obtener una descripción más detallada, consulte el capítulo 4 Asistente Crear o editar método.

Iconos

21

Icono: Cambiar usuario actual

Si la administración de usuarios está activa (consulte el capítulo 9.4 Administración de usuarios (magellan Tracker) y el capítulo 9.5 Administración de usuarios (magellan Standard)) haga clic en este icono para cerrar la sesión del usuario actual y volver a entrar con otro usuario.

Con **magellan** Standard la administración de usuarios es opcional. Con **magellan** Tracker la administración de usuarios es obligatoria.

🚽 Icono: Otras opciones

haga clic en el icono Otras opciones para seleccionar las opciones siguientes:

	Control del instrumento		Opciones
	Procesamiento de archivos	22	Administración de usuarios
m	Acerca de magellan		

 El Control del instrumento ofrece un rápido acceso a diferentes funciones del instrumento, así como a las opciones de configuración y servicios. Consulte el capítulo 3 Control y configuración del instrumento.

 El Procesamiento de archivos se utiliza para mover archivos a una ubicación de archivo, convertir archivos a o desde otra versión de magellan, o para importar datos sin procesar de un archivo ASCII. Consulte el capítulo 9.2 Procesamiento de archivos.



- Opciones se utiliza para personalizar ciertos ajustes predeterminados como el directorio de los archivos generados, las opciones de portapapeles y copia Excel, la vista de placa y los ajustes del asistente otras opciones, inicio, idioma y contraseña. Consulte el capítulo 9.3 Opciones.
- La Administración de usuarios se utiliza para añadir o eliminar usuarios y para establecer o modificar derechos de usuario. Consulte el capítulo 9.4 Administración de usuarios (magellan Tracker), y respectivamente el capítulo 9.5 Administración de usuarios (magellan Standard).
- Acerca de magellan ofrece información de licencias y detalles sobre la versión y los componentes actualmente instalados. El registro se puede solicitar mediante el Asistente para registrar (consulte el capítulo 2.4.3 Registrar magellan).

Cierre la ventana Otras opciones para regresar a la lista de asistentes.

Icono: Control de temperatura

Utilice este icono para establecer la temperatura objetivo del instrumento conectado.

Para obtener una descripción más detallada, consulte el capítulo 3.1.2 Control de temperatura....



Icono: Mover la placa

Utilice este icono para mover el portaplacas dentro o fuera del instrumento.

2.2 Componentes y términos – Lógica básica de magellan

2.2.1 Tipos de archivo usados con magellan

Se utilizan los tipos de archivo siguientes con magellan.

Por defecto, todos los tipos de archivo asociados con **magellan** se guardan en subdirectorios en el directorio adecuado:

...\All Users\Documents\Tecan\

Los subdirectorios se muestran en la siguiente tabla:

Tipo de archivo	Extensión de archivo	Directorio
Zona de trabajo	.wsp	\magellan\wsp
Método	.mth	\magellan\mth
Lista de IDs de muestras	.smp	\magellan\smp
Archivos de exportación	.asc	\magellan\asc
Curva estándar	.std	\magellan\wsp
Definición de placa	.pdf / pdfx	\Reader\pdf
		\pdfx

Para obtener una descripción detallada de los métodos, listas de IDs de muestras, zonas de trabajo y curvas estándar, consulte los capítulos 4 Asistente Crear o editar método, 5.2.3 Importar lista de IDs de muestras y 7 Asistente Evaluar resultados.



Aviso Los instrumentos INFINITE F50 utilizan el formato de archivo .pdfx, mientras que los instrumentos SUNRISE utilizan archivos .pdf.

TECAN.



Aviso Los menús y barras de herramientas disponibles cambian en función del tipo de archivo abierto.

2.2.2 Procesamiento de carpetas

En **magellan** es posible crear una estructura de carpetas específica de usuario para guardar archivos (métodos, zonas de trabajo, curvas estándar y listas de ID de muestra) en cualquier carpeta del Explorador de Windows. Las carpetas nuevas se pueden crear haciendo clic con el botón secundario del ratón y seleccionando **Nueva carpeta** en el menú contextual. Los archivos y carpetas se pueden mover/cortar/copiar de la misma manera que con el Explorador de Windows.

Símbolos de tipo de archivo

magellan utiliza los siguientes símbolos para métodos, zonas de trabajo, curvas estándar y listas de ID de muestra en la ventana de vista de archivo:

0	Método
	Curva estándar
6	Zona de trabajo
<u></u>	Método protegido por contraseña
P	Lista de IDs de muestras

magellan Standard

Los directorios predeterminados para guardar los archivos creados nuevos se pueden establecer mediante: página principal de Lista de asistentes \rightarrow botón **Otras opciones** \rightarrow botón **Opciones** \rightarrow pestaña **Directorios**.

Los usuarios pueden crear carpetas nuevas en cualquier carpeta del Explorador de Windows durante el proceso de guardado.

En la ventana **Guardar**, la carpeta, especificada por el directorio predeterminado del archivo, se abre automáticamente cuando se crear un archivo nuevo. Si se modifica un archivo existente, se abre el directorio de la ubicación actual. Sin embargo, es posible guardar archivos en cualquier carpeta del Explorador de Windows o en una carpeta nueva.



magellan Tracker

Los usuarios de **magellan** con derechos de administración pueden establecer los directorios predeterminados para guardar nuevos archivos creados mediante: página principal de **Lista de asistentes** → botón **Otras opciones** → botón **Opciones** → **Directorios**. Estos directorios predeterminados son válidos para todos los usuarios.

Los usuarios pueden crear nuevas subcarpetas solo si tienen privilegios de administrador de Windows y solo en la ruta especificada como predeterminada durante el proceso de guardado.

En la ventana **Guardar**, la carpeta, especificada por el directorio predeterminado del archivo, se abre automáticamente cuando se crear un archivo nuevo. Si se modifica un archivo existente, se abre el directorio de la ubicación actual. Sin embargo, solo se pueden guardar archivos en la carpeta predeterminada o en una subcarpeta nueva o existente dentro de esta carpeta.

2.2.3 Elementos estándar

Cada asistente muestra secuencialmente un número de ventanas en las que se ofrece toda la información necesaria, ajustes y posibilidades de entrada de datos.

Elementos estándar de los asistentes de Magellan

Para navegar de una ventana a la otra existen varios botones en la parte inferior de la página del asistente:

Botón Atrás	El botón Atrás se utiliza para regresar a la ventana previa dentro de un asistente.
Botón Siguiente	El botón Siguiente se utiliza para navegar a la ventana siguiente dentro de un asistente.
Botón Finalizar	El botón Finalizar se utiliza para ir a la ventana de guardar del asistente.
Botón Guardar	El botón Guardar solo se encuentra en la última ventana del asistente y sustituye al botón Siguiente . Se utiliza para cerrar el asistente y guardar todos los cambios o para iniciar un proceso.
Botón Cancelar	El botón Cancelar se utiliza para cerrar un asistente sin guardar los cambios en los ajustes o documentos.
Botón Ayuda	El botón Ayuda abre la ventana de ayuda.

Elementos estándar de Microsoft Windows

Botón Aceptar	Este botón confirma los ajustes, aplica y guarda los cambios consecuentemente y cierra el cuadro de diálogo.
Botón Cancelar	Este botón cierra el cuadro de diálogo sin guardar ningún cambio en los ajustes o documentos.
Botón Ayuda	Haga clic en el botón Ayuda para abrir la ayuda en línea de magellan.

Información de barra de estado

La barra de estado muestra la información siguiente:

- Información del comando actual
- Nombre del usuario con la sesión actualmente abierta.
- Nombre del instrumento conectado. Por ejemplo: Sunrise



- Método: modo de medición y unidad. Por ejemplo: Absorbencia [DO]
- Zona de trabajo: fecha y hora de la medición. Por ejemplo: 27.11.2002 14:13:03
 - Número de pocillos seleccionados. Por ejemplo: 3 pocillo(s) seleccionado(s)
- Información de estado del teclado: actividad de las teclas de alternar del teclado: MAY (bloqueo de mayúsculas), NUM (bloqueo del teclado numérico), DESP (bloqueo de desplazamiento)
- Icono de estado de conexión de instrumento

2.2.4 El botón de Ayuda

Haga clic en el botón **Ayuda** o pulse 'F1' para abrir la ayuda en línea de **magellan**.

2.2.5 El cuadro de diálogo de bienvenida

Cada asistente empieza con un cuadro de diálogo de **Bienvenida**, que ofrece una breve descripción introductoria del procedimiento a realizar.

Desmarque la casilla de verificación **Mostrar página de bienvenida** para suprimir las páginas de bienvenida al iniciar los asistentes en el futuro.

2.2.6 Lista de accesos directos

MAYÚS+B	Botón Atrás
MAYÚS+N	Botón Siguiente
ESC	Botón Cancelar
INTRO	Botón Siguiente o Finalizar, si la ventana está activa
F1	Menú Ayuda
CTRL+C o CTRL+INSERTAR	Copiar
CTRL+V o MAYÚS+INSERTAR	Pegar
CTRL+X	Cortar
CTRL+Y	Rehacer
CTRL+Z	Deshacer
SUPR	Eliminar el contenido del pocillo activo (editar ID de muestra, editar fórmula)
CTRL+MAYÚS	Mostrar la fórmula del pocillo seleccionado al visualizar el resultado de la transformación (asistente Evaluar resultados)



2.3 Iniciar magellan

2.3.1 Iniciar la versión estándar

Realice los pasos siguientes para iniciar **magellan**:

- 1. Asegúrese de que se ha conectado un instrumento o seleccione un instrumento simulado.
- 2. En el menú **Inicio** de Windows, seleccione el grupo de programas **Tecan** y haga clic en el icono magellan.
- 3. magellan se inicia.



Aviso magellan se puede ejecutar con un instrumento conectado o en modo de demostración (simulando un instrumento). Si desea conectar un instrumento (consulte el capítulo 2.4 Conectar un instrumento), encienda el instrumento antes de iniciar magellan.

Aviso Antes de empezar a trabajar con magellan, se recomienda leer el documento Release Notes for magellan por cualquier mal funcionamiento de la aplicación.

2.3.2 Iniciar la versión Tracker



Aviso
Al utilizar magellan Tracker, el sistema de administración de usuarios
debe personalizarse mediante un administrador designado, responsable
de la configuración de las cuentas de usuario y de la asignación de
derechos de usuario. El administrador siempre debe haber sido formado
por Tecan o una organización autorizada por Tecan.

Iniciar sesión en el sistema por primera vez con magellan Tracker

Al iniciar **magellan Tracker** por primera vez, aparece un cuadro de diálogo, informando al usuario de que primero hay que crear un **Administrador de usuarios**. Haga clic en **OK** y aparecerá el cuadro de diálogo **Crear administrador**.

Rellene los campos de texto y haga clic en **Aceptar** para guardar los ajustes. Como mínimo debe crearse un administrador de usuarios.



PRECAUCIÓN SI SE OLVIDA LA CONTRASEÑA DEL ADMINISTRADOR DE USUARIOS, NO SE PUEDEN MODIFICAR LOS AJUSTES DE LA ADMINISTRACIÓN DE USUARIOS Y MAGELLAN DEBE REINSTALARSE POR COMPLETO.

RECOMENDAMOS CREAR COMO MÍNIMO DOS ADMINISTRADORES DE USUARIOS.



Para más definiciones de usuario de la administración de usuarios (p. ej. derechos de usuario), consulte el capítulo 9.4 Administración de usuarios (magellan Tracker) y el capítulo 9.5 Administración de usuarios (magellan Standard).

Contraseña

Contraseña inicial (solo magellan Tracker)

Cuando un usuario entra en el sistema por primera vez, hay que cambiar la contraseña asignada por el administrador.

Después de introducir el **Nombre de usuario** y la **Contraseña** y hacer clic en **Aceptar**, se mostrará el mensaje "Su contraseña inicial solo es válida para una ocasión. ¡Debe cambiar la contraseña!". Haga clic en **Aceptar**; la ventana **Cambiar contraseña** obliga al usuario a introducir una contraseña nueva.

Caducidad de la contraseña (solo magellan Tracker)

Cuando caduca la contraseña, se pide al usuario que introduzca una contraseña nueva.

Consulte el capítulo 9.4.4 Opciones de Administración de usuarios.

Inicio de sesión

Una vez que se haya activado la administración de usuarios, aparecerá el cuadro de diálogo **Inicio de sesión** cada vez que se inicie **magellan**.

El cuadro de diálogo Inicio de sesión contiene los siguientes elementos:

Cuadro de texto Nombre de usuario	Introduzca el ID de usuario
Cuadro de texto Contraseña	Introduzca la Contraseña

Haga clic en el botón IR para entrar o en el botón Cancelar para cerrar magellan.

Aplicación bloqueada

Si la aplicación no se utiliza durante un tiempo máximo especificado (definido por el usuario únicamente en **magellan** Tracker), entonces se bloqueará. Se deberá escribir la contraseña para desbloquear la aplicación.



2.4 Conectar un instrumento

2.4.1 Conectarse a instrumentos SUNRISE

Conecte el instrumento al ordenador según se describe en las Instrucciones de uso del instrumento y encienda el instrumento antes de iniciar **magellan**.

Primer inicio de magellan

La primera vez que se inicia **magellan** se mostrará el cuadro de diálogo **Instalar puerto**:

Instrumento Seleccionar:	SUNRISE	Aceptar Cancelar
Puerto		Ayuda
 Seleccionar: Buscar cualquiera Modo de demostr 	Buscar cualquiera ~	
nstrumento	Seleccione el instrumento conectado de la o seleccione Buscar cualquiera .	ı lista desplegat
uerto	Seleccione el Puerto COM adecuado de la o seleccione Buscar cualquiera y haga c	a lista desplega lic en Aceptar .

Cambiar instrumento

	_	_	_	
- 6	2	-	_	ł
	۱.,	- 11	14	I
	Ŀ			

Haga clic en el icono **Otras opciones** en la lista de asistentes y en **Control del instrumento** en la lista de otras opciones. Seleccione **Cambiar instrumento** en **Configuración & Servicios** para conectar un instrumento o para cambiar el instrumento actualmente conectado.

Modo demostración

Si no hay **ningún instrumento** conectado, seleccione el instrumento a simular de la lista desplegable en el cuadro **Instrumento** y seleccione **Modo demostración** en el cuadro de **Puerto** y haga clic en **Aceptar**. Ahora **magellan** está en modo demostración.

Instrumento			Aceptar
Seleccionar:	SUNRISE	~	
O Buscar cualquiera			Cancelar
Puerto			Ayuda
O Seleccionar:	Buscar cualquiera	\sim	
O Buscar cualquiera		51	
Modo de demostr	ación		

Modo demostración permitido (en el cuadro **Opciones**) está activado de modo predeterminado.

En el modo de demostración es posible realizar todas las funciones excepto ejecutar una medición real.

2.4.2 Conectarse a instrumentos INFINITE F50

Primer inicio de magellan

La primera vez que se inicia **magellan** se mostrará el cuadro de diálogo **Instalar puerto**:

nstrumento			Aceptar
Seleccionar:	infinite F50	~	
) Buscar cualquiera	а.		Cancelar
uerto			Ayuda
) Seleccionar:	Buscar cualquiera	.*	

Seleccione **INFINITE F50** en el cuadro de grupo **Instrumento** del cuadro de diálogo **Instalar puerto**. Haga clic en **Aceptar** para mostrar el cuadro de diálogo **Conectar al instrumento**, seleccione el instrumento y haga clic en **Aceptar**.

Conectar a:			
Nombre del instrumento	Тіро	Alias	Puerto
infinite F50	READER	160 100 36 20	USB0
Conexión adicional a:			
Nombre del instrumento			Puerto
] Mostrar los instrum <mark>e</mark> ntos sir	nulados		

Seleccione el instrumento y haga clic en Aceptar.



Modo Demo

Si **no hay ningún instrumento** conectado, seleccione en la lista desplegable el instrumento que se va a simular en el cuadro de grupo **Instrumento** del cuadro de diálogo **Instalar puerto** y haga clic en **Aceptar**.

onectar al instrumento			
Conectar a:			
Nombre del instrumento	Tipo	Alias	Puerto
infinite F50	Reader	Simulation	SUSIM:
Conevién adicional au			
Nombre del instrumento			Puerto
ConnectSimulator			Connec
7 Montrar los instrumentos sin	nulados		
		1.30_1111163	,
Acer	otar		Cancelar

Para el modo de demostración, seleccione la casilla de verificación **Mostrar instrumentos simulados** y el instrumento preferido en la lista desplegable y haga clic en Aceptar.

2.4.3 Registrar magellan



Aviso

magellan se puede usar sin registrar durante 30 días de trabajo. Si magellan no se registra tras este período, las opciones de guardar e imprimir se desactivan.

El registro de **magellan** se realiza con el asistente **Registro** de **magellan**. Si **magellan** no se ha registrado, este asistente se inicia automáticamente cada vez que se abre el programa.

El asistente **Registro** de **magellan** también se puede abrir mediante el icono **Otras opciones** en la página de inicio de la lista de asistentes. (**Otras opciones** \rightarrow **Acerca de magellan** \rightarrow **Registro**).

Para registrar **magellan**, se necesita un número de serie de Magellan, el número de identificación de unidad de hardware (HUID) y un número de licencia.

Existen dos posibilidades para registrar magellan:

1. El software se ha comprado. Junto con el software se suministra el número de serie.

Después de introducir el número de serie en el cuadro de diálogo de registro de **magellan**, rellene el formulario de registro para iniciar la solicitud del número de licencia. El formulario de registro rellenado se envía junto con el HUID individual



a Tecan Austria. El software genera el número HUID y lo introduce automáticamente en el formulario de registro. Está relacionado con el número de unidad de sistema ofrecido por Windows.

Tras recibir el número de licencia debe iniciar el asistente de registro de nuevo e introducir el número. El asistente **Registro** de **magellan** confirma el número de licencia y resume la información del usuario. Haga clic en el botón **Finalizar** para completar el procedimiento de registro. Entonces las funciones de **magellan** estarán totalmente disponibles para el usuario.

 El software se ha suministrado como versión de demostración junto con el instrumento. Posteriormente es posible un registro (no se envía un número de serie).

Para adquirir magellan y obtener un número de serie, seleccione la opción **Haga un pedido de Magellan para obtener una licencia válida** y siga con el procedimiento de registro según se ha indicado antes.



Aviso Es de suma importancia que la persona que registra el software también tenga derechos de administrador en el sistema operativo del ordenador.

2.4.4 Asistente para registrar

El asistente para registrar se inicia con el cuadro de diálogo de **Bienvenida**, que contiene una breve descripción del asistente. Haga clic en **Siguiente** y se mostrará el cuadro de diálogo **Número de serie**.

Números de serie y de licencia

Para solicitar un número de serie para magellan

Para adquirir **magellan** y recibir un número de serie, seleccione la opción **Haga un pedido de Magellan para obtener una licencia válida**. Se abrirá la ventana **Formulario de registro** (consulte la sección Formulario de registro a continuación).

Registrar Magellan	×
Escriba su número de serie Magellan:	
O Haga un pedido de Magellan para obtener una licencia válida.	
C Dank Navit	<u> </u>



Si ya ha adquirido magellan

Seleccione **Escriba su número de serie magellan**. Anote su número de serie incluido en el paquete de **magellan**.

Después de introducir el número de serie, haga clic en **Siguiente** y se mostrará el cuadro de diálogo **Número de licencia**.

Escriba su número de licencia Magellan:	
Solicita un número de licencia Magellan porque no la recibió.	

Los usuarios que instalen el software por primera vez no contarán aún con un número de licencia y deberán seleccionar **Solicita un número de licencia Magellan porque no lo recibió**. Haga clic en **Siguiente** y se abrirá la ventana **Formulario de registro**.

Formulario de registro

Complete el formulario de registro. Los campos marcados con un asterisco son obligatorios.

Nombre completo:	Test User				
Empresa / Institución:	Tecan Austria				
Dirección postal 1:	Untersbergstrasse	1a			•
Dirección postal 2:					
Código postal, Población:	5082 Groedig	* País:	Austria	~	•
Teléfono:	I	Fax:			
Correo electrónico:	test.user@tecan.co	om		2	
Área de trabajo (Ciencias Diagnóstico cl <i>í</i> nico /Otra:	de la vida /			~	
Instrumentos Tecan utilizados:	SUNRISE			Añadir	
Detection Modes	Absorbance			Añadir	
	Quiero material ir	nformativo de Teca	an.		
		12 12	87	-214	



Haga clic en **Siguiente** y se abrirá la ventana de Contrato de licencia. Léalo detenidamente y haga clic en el botón **Acepto**.

Haga clic en Siguiente y se mostrará la información de registro.

USER ADDRESS: Test Unter	User / Tecan Austria sbergstrasse la	^
5082 APPLICATION: Magel WORKSTATION: 9408G	Groedig / Austria lan, 7, 2, 1, 5 -3U335FB	
<		>
mulario de correo electrónico	Imprimir formulario	Ver formulario.

Haga clic en **Formulario de correo electrónico...** para enviar la información a través de su programa de correo electrónico predeterminado de su ordenador o bien pulse **Imprimir formulario...** para imprimirlo y enviarlo por correo o fax a Tecan. Haga clic en **Ver formulario...** para visualizar el formulario de registro en el Bloc de notas o en WordPad. Dentro de las siguientes 24 horas recibirá el número de licencia.

Completar la licencia de magellan

Después recibir el número de licencia de Tecan, repita los pasos anteriores e inserte el número de licencia. Haga clic en **Siguiente**. El asistente de registro entonces confirmará el número de licencia y hará un resumen de la información del usuario.

Haga clic en **Finalizar**. Se cerrará el asistente **Registro de magellan** y las funciones de **magellan** quedarán a plena disposición del usuario.



3. Control y configuración del instrumento

Haga clic en el botón **Control del instrumento** en el cuadro de diálogo **Lista de asistentes** y se mostrará el cuadro **Control del instrumento**.

En función del instrumento conectado se activan o desactivan diferentes opciones de configuración y servicios del instrumento.

3.1 Opciones del instrumento

3.1.1 Movimientos...

Esto abre el cuadro de diálogo **Movimientos**, en el cual es posible controlar los movimientos del portaplacas y del placa de filtro.

3.1.2 Control de temperatura...

Solo disponible para instrumentos equipados con control de temperatura.

Esta opción permite al usuario establecer la temperatura dentro del lector.

Este cuadro también está disponible a través del cuadro de diálogo **Iniciar medición** antes de iniciar una medición. (Consulte el capítulo 6.5 Iniciar medición con un método predefinido o favorito).

Control de tem	peratura	
Temperatura ac Temperatura ob Cor	tual Temperatura actual: 32.2 °C Recargar jetivo atrol de temperatura activo Temperatura objetivo: 37 °C Aplicar	Aceptar Cancelar Ayuda Introducir placa
Temperatura actual	Se muestra la temperatura ac correspondiente. Haga clic er actualizar la temperatura actu	ctual en el campo n el botón Recargar para ual.
Temperatura	Para iniciar o detener el conti	rol de temperatura, marque o

objetivodesmarque la casilla de verificación e introduzca la
temperatura objetivo. Haga clic en Aplicar
para enviar la temperatura al instrumento y luego haga clic
en Aceptar para cerrar el cuadro de diálogo.

3.2 Opciones de Configuración y Servicios

3.2.1 Cambiar instrumento...

En el cuadro **Configuración y Servicios**, haga clic en **Cambiar instrumento...** para abrir el cuadro de diálogo **Instalar puerto**.

Esta opción permite conectar **magellan** a un instrumento. Para obtener más detalles, consulte el capítulo 2.4 Conectar un instrumento.



Aviso

Si conecta otro instrumento al ordenador o si modifica los parámetros de la interfaz, siempre deberá seleccionar esta opción de menú. Los ajustes se utilizan automáticamente la próxima vez que se inicia el software.

3.2.2 Definir placas de filtro...

Permite definir placas de filtro estándar y definidas por el usuario.

3.3 Archivos de registro

Durante el funcionamiento de **magellan** se crean archivos de registro. La comunicación entre el software y el instrumento y la comunicación entre los componentes del software se almacena en estos archivos de registro.

Los archivos se encuentran en los directorios siguientes:

- Archivos de registro de magellan (comunicación entre los componentes individuales de magellan): Windows 10: C:\Users\Public\Documents\Tecan\LogFiles\
- Conjunto de archivos de registro necesarios
 Los archivos de registro se pueden guardar como archivo zip seleccionando
 el botón Guardar archivos de registro... en el cuadro de diálogo
 Procesamiento de archivos (Otras opciones Procesamiento de
 archivos). Entonces podrá escribir un nombre para el archivo zip y guardarlo
 en un directorio definido. En caso de cualquier error de estado o de medición
 de Magellan, este archivo también contiene datos, estados (p. ej.
 desbordamiento, lámpara tenue) o errores de cálculo, y se puede enviar
 fácilmente a su departamento de soporte local para solicitar soporte. Para
 obtener información adicional, consulte el capítulo 9.2 Procesamiento de
 archivos Guardar archivos de registro.
- Archivos de registro de INFINITE F50 (comunicación entre magellan e instrumentos INFINITE F50): Windows 10: C:\Users\Public\Documents\Tecan\LogFiles\Magellan\V x.y\Instrument Serial Number
- Archivos de registro SUNRISE (RdrOle.log; comunicación entre magellan e instrumentos SUNRISE) de modo predeterminado: Windows 10: C:\Users\CurrentUser\AppData\Local\Temp



4. Asistente Crear o editar método

4.1 Introducción

El asistente Crear o editar método se utiliza para

- Crear o editar métodos,
- Fijar parámetros de medición y evaluación,
- Definir el diseño de placa,
- Seleccionar el formato del informe impreso y
- Fijar los parámetros del procesamiento automatizado de datos.

Resumen del flujo de trabajo

Haga clic en **Continuar** en la página de bienvenida del asistente **Crear o editar método**. En la ventana siguiente, seleccione

- Crear nuevo si desea definir un método nuevo o
- Editar si desea modificar un método existente.

En la ventana **Definir parámetros de medición** se pueden fijar parámetros de medición. Haga clic en **Definir evaluación** para definir el diseño de placa, parámetros de evaluación, parámetros de impresión y procesamiento de datos automatizado. Al final del asistente el método nuevo o modificado se guarda como archivo .mth.



Aviso Para obtener una descripción paso a paso acerca de cómo crear un método, consulte 12 Ejemplo de aplicación.

Página de selección de archivo

En la lista de asistentes, haga clic en el botón **Crear o editar método**. Haga clic en **Siguiente** en la página de bienvenida del asistente **Crear o editar método** y aparecerá la página de **Selección de archivo** con los elementos siguientes:

Botón Crear nuevo	Seleccione el botón Crear nuevo para crear un método nuevo.
Botón Editar	Debe seleccionar el botón Editar para editar un método existente.
Cuadro combinado Mostrar	En el cuadro combinado Mostrar se puede definir un filtro de archivo para que en la lista solo aparezcan los métodos que cumplen los criterios seleccionados:
	Todos los archivos
	Archivos de este instrumento
	 Mis archivos: esta opción solo estará disponible si está habilitada la administración de usuarios (siempre está habilitada en magellan Tracker).
	Archivos firmados: solo disponible para magellan Tracker
	 Archivos de ejemplo: solo estarán disponibles si se han instalado.

Lista Nombre de	Seleccione el método a editar desde la lista Nombre de archivo .
archivo	Un campo de Notas junto a cada nombre de archivo contiene, si se ha escrito, una breve descripción del método.
	Se muestran todos los métodos disponibles en el directorio de métodos estándar (consulte el capítulo 9.3 Opciones).
Botón Presentación preliminar	Haga clic en el botón Presentación preliminar para abrir el cuadro de diálogo Presentación preliminar , donde se muestra una vista previa de los ajustes del método seleccionado y se puede iniciar una impresión.

TECAN.



Aviso

Si tiene un instrumento SUNRISE conectado, pero selecciona un método de INFINITE F50, se mostrará el cuadro de diálogo de Incongruencia de instrumento:

Conecte el instrumento adecuado

4.2 Definir los parámetros de medición

En la ventana **Parámetros de medición** es posible fijar todos los parámetros necesarios para la medición, incluyendo el modo de medición, longitudes de onda, modo de lectura, temperatura, etc. en función del tipo de instrumento conectado.

Si tiene conectado un instrumento INFINITE F50, para conocer más detalles acerca de cómo definir los parámetros de medición, consulte las Instrucciones de uso de i-control.



Aviso Los parámetros disponibles a definir dependen del instrumento conectado.



4.2.1 Tipos de mediciones - SUNRISE

Haga clic en **Haga su selección** y se mostrará el cuadro de diálogo **Tipo de medición**:

Crear o editar métod	0		×
	Tipo de l	medición Medición de punto final Medición cinética Medición cinética de pocillo Medición de múltiples etiquetas Parámetros de medición	
Ayuda			
Cancelar	<<< Atrás		
		© 2016 Tecan	

El tipo de medición se debe seleccionar en una lista de botones de opciones.

La ventana **Tipo de medición** contiene los siguientes elementos (en función del instrumento conectado):

Cuadro de grupo Tipo de medición	•	Seleccione Medición de punto final para realizar una medición simple.
	•	Seleccione Medición cinética para realizar mediciones cinéticas por placa con un intervalo de tiempo especificado.
	•	Seleccione Medición de etiquetas múltiples para realizar mediciones de etiquetas múltiples con distintos parámetros de medición.

Haga clic en **Parámetros de medición...** para abrir el cuadro de diálogo Parámetros de medición.

4.2.2 Parámetros de medición - SUNRISE

Seleccione el **Tipo de medición** requerida y haga clic en **Parámetros de medición...** para abrir el cuadro de diálogo Parámetro de medición para ajustar

- Códigos de barras
- Longitudes de onda
- Parámetros de medición (modo lectura)
- Cinéticas
- Temperatura y
- Agitación.

Haga clic en el botón **Aceptar** para guardar los cambios o haga clic en el botón **Cancelar** para rechazarlos.

Según el instrumento conectado y el tipo de medición seleccionado, varían el cuadro de diálogo **Parámetros de medición** y las pestañas respectivas disponibles.



Cuando se conecte a un instrumento Infinite, consulte las Instrucciones de uso para i-control.

Medición de etiquetas múltiples

Si se selecciona el tipo de medición de etiquetas múltiples, se muestra el cuadro de diálogo **Medición de etiquetas múltiples**.

Para crear una lista de ventanas de parámetros deberá hacer clic en el botón **Nuevo**. Se mostrará el cuadro de diálogo **Parámetros de medición** para definir un nuevo conjunto de parámetros de medición. En el cuadro de diálogo de etiquetas múltiples se deben crear al menos dos conjuntos de parámetros de medición.



Aviso

Si al ejecutar el método se selecciona Sacar placa después de la medición, magellan mostrará un cuadro de mensaje que permite modificar los líquidos en la placa y continuar con la medición. De lo contrario, la medición se realiza sin ninguna interrupción.

El parámetro de medición **Comentarios** se puede usar para etiquetar el conjunto de parámetros de medición de la lista.

El cuadro de diálogo Etiquetas múltiples contiene los siguientes elementos:

I DEF-1 Modo de medición: Absorbencia;Lor 2 DEF-2 Modo de medición: Absorbencia;Lor 3 DEF-3 Modo de medición: Absorbencia;Lor Borrar Cinética Número de ciclos: 2 Intervalo: 000:00:00 hh:mm:ss Usar intervalo mínimo	Núm.	Etiqueta	Parámetros de medición	Nuevo
2 DEF-2 Modo de medición: Absorbencia;Lor 3 DEF-3 Modo de medición: Absorbencia;Lor Borrar Cinética Número de cidos: 2 Intervalo: 000:00:00 hh:mm:ss Usar intervalo mínimo	1	DEF-1	Modo de medición: Absorbencia;Lo	Propiedades
Sorrar Modo de medicion: Absorbencia;Lor Borrar Cinética Número de cidos: 2 Intervalo: 000:00:00 hh:mm:ss Usar intervalo mínimo	2	DEF-2	Modo de medición: Absorbencia;Lo	r -
< Cinética Número de cidos: Intervalo: Usar intervalo mínimo				borran
Cinética Número de cidos: 2 Intervalo: 000:00:00 hh:mm:ss Usar intervalo mínimo				
Cinética Número de ciclos: 2 Intervalo: 000:00:00 hh:mm:ss Usar intervalo mínimo				
Número de ciclos: 2 Intervalo: 000:00:00 hh:mm:ss Usar intervalo mínimo	<		>	
Intervalo: 000:00:00 hh:mm:ss	< Cinétio	ca	>	
Usar intervalo mínimo	< Cinétio Núr	ca nero de ciclos:	2	
	< Cinéti Núr Inte	ca nero de cidos: ervalo:	2 000:00:00 hh:mm:ss	

Lista Etiquetas En la lista de **Etiquetas múltiples** se enumeran línea a línea las definiciones de parámetros de medición existentes. En la columna **Núm.** un número guía cuenta las definiciones de parámetros de medición existentes y se muestra un pequeño icono. La columna **Etiqueta** muestra un nombre personalizable (en **Parámetros de medición – Pestaña General – Comentario a esta medición**). La columna **Parámetros de medición** enumera un resumen de los parámetros de medición seleccionados.


Botón Nuevo	Haga clic en el botón Nuevo y podrá definir los parámetros de medición de cada nueva medida. Se mostrará el cuadro de diálogo Parámetros de medición .
	Defina nombres para cada etiqueta. En el cuadro de diálogo Parámetros de medición se utiliza como etiqueta el nombre escrito en el campo de comentario de la pestaña General .
	Haga clic en el botón Nuevo y podrá definir los parámetros de medición de cada nueva medida. Se mostrará el cuadro de diálogo Parámetros de medición .
	Defina nombres para cada etiqueta. En el cuadro de diálogo Parámetros de medición se utiliza como etiqueta el nombre escrito en el campo de comentario de la pestaña General .
	Para definir una nueva medición deberá introducir todos los parámetros.
Botón Propiedades	Haga clic en el botón de Propiedades y se mostrarán y se podrán editar todos los parámetros de medición de la etiqueta resaltada en la lista de parámetros de medición.
Botón Eliminar	Haga clic en Borrar o pulse la tecla Supr para borrar los parámetros de medición resaltados del campo de etiquetas múltiples.
Número de ciclos	Define el número de ciclos cinéticos para una medición cinética de etiquetas múltiples.
Intervalo	Define el intervalo cinético o la pausa entre etiquetas (si solo se define un ciclo).

4.2.3 Parámetros de medición - INFINITE F50

Definición de mediciones de punto final

El siguiente ejemplo describe una **Medición de punto final de absorbencia** en todos los pocillos de una microplaca de 96 pocillos.

- 1. Seleccione una microplaca de 96 pocillos en la lista desplegable **Definición de placa**.
- 2. De modo predeterminado, se selecciona la medición de todos los pocillos de la microplaca de 96 pocillos.
- 3. Escriba la medición que desee y las longitudes de onda de referencia.

	2
6	
	3
Ebqueta	
Nompre: Laber	
alb	Bouch Nontre Lodi v



Definición de mediciones de etiquetas múltiples

El siguiente ejemplo describe una **Medición de etiquetas múltiples de absorbencia** en un rango definido de una microplaca de 96 pocillos (A1:E7). Se medirán tres etiquetas de absorbencia.

- 1. Seleccione una microplaca de 96 pocillos en la lista desplegable **Definición** de placa.
- 2. De modo predeterminado, se selecciona la medición de todos los pocillos de la microplaca de 96 pocillos.
- 3. Escriba la longitud de onda de medición que desee.
- 4. Inserte dos elementos más de **Absorbencia** y escriba las longitudes de onda de medición.

🐐 🔻 Sección de placa		2
	ale	
🛔 🖛 Absorbencia		3
- Longitud de anda Medición: 405 nm ✓ ☐ Referencia: 405 nm ✓	Etiqueta Nombre: Etiqueta 1 v	
🛔 🔻 Absorbencia		4
Longitud de onda Medición: 492 nm ~ Referencia: 405 nm ~	Ekgeta Nombre Ekgeta2 v	
🕴 🖛 Absorbencia		5
- Longitud de onda Medición: 620 nm ✓ □ Referencia: 405 nm ✓	Eliqueta Nombre: [Eliqueta3 v	

Definición de mediciones cinéticas

El siguiente ejemplo describe una medición cinética de una microplaca de 96 pocillos.

- 1. Seleccione una microplaca de 96 pocillos en la lista desplegable **Definición** de placa.
- 2. Inserte un elemento de programa **Ciclo cinético** entre la sección de placa y el elemento de absorbencia.
- 3. Ciclos/Número de ciclos: 50
- Intervalo cinético (intervalo entre mediciones): seleccione Utilizar intervalo cinético y escriba: 2 minutos 30 segundos.
- 5. Escriba la longitud de onda de medición que desee para definir el elemento **Absorbencia**.

🔹 🔻 Sección de placa		2
123441471180000 000000000000000000000000000000000000	Detales_	
😡 🔻 Ciclo cinético		3
Cidos	Intervalo cinético	
Número de ciclos: 50 ÷	Utilizar intervalo cinético:	
O Duración	(0) 022:30	
Absorbencia		4
Longitud de onda	Etiqueta	
Medición: 492 nm	Nombre: Etiqueta1 V	
Referencia: 405 rm		



4.3 Definir evaluación

4.3.1 Ventana de visión global de Crear o editar método

En la ventana **Definir evaluación** el usuario define el diseño de placas, las transformaciones y los cálculos, selecciona el formato del informe impreso y establece los parámetros de procesamiento automatizado de fechas.



Barra de herramientas

Encima de la ventana hay una **Barra de herramientas** con las funciones más comunes en función de las opciones actualmente seleccionadas:

Botón Deshacer	Haga clic en Deshacer para anular una acción previa.
Botón Rehacer	Haga clic en Rehacer para repetir una acción que se había anulado.
Botón Seleccionar todas las placas sin usar	Solo en el modo Diseño de placas: Se seleccionan todos los pocillos sin usar de la placa.
Botón Zoom al 10%	Fija la vista del diseño de placas al 10%.
Botón Zoom al 100%	Fija la vista del diseño de placas al 100%.
Botón Modo zoom	Utilice el botón Modo zoom para ampliar el área marcada. Cuando se activa, el usuario puede seleccionar un área para ampliar haciendo clic y arrastrando un marco sobre el área de diseño deseada. Haga clic con el botón secundario del ratón para aleiar la vista al 100%.



Ventana de diseño de placas

En el área central de la ventana, la **Ventana de diseño de placas** muestra un diseño esquemático de una microplaca.

Al crear un método nuevo, seleccione los pocillos para la medición. Un clic de ratón selecciona un pocillo individual. Los pocillos no seleccionados permanecen en gris.

Para el diseño definido, los identificadores, transformaciones y fórmulas seleccionadas en la **Barra de control** aparecen en los pocillos correspondientes (consulte **Barra de control de Crear o editar método** más adelante).

Diseño de la placa: menú contextual

Haciendo clic con el botón secundario del ratón en el diseño de placas aparece un menú **contextual** para los pocillos marcados. Están disponibles los siguientes comandos:

Menú	Descripción
Resumen	En todas las vistas de placa. Se muestra el cuadro de diálogo Resumen . Consulte en la sección 7.4.18 Menú contextual de un pocillo/ Cuadro de diálogo Resumen información adicional sobre el cuadro de diálogo Resumen Esta opción está disponible si se ha asignado un diseño al pocillo seleccionado.
Selección de relleno	Solo en el modo Diseño de placas. Si se ha marcado un pocillo o un área de la placa, puede rellenarse con los identificadores correspondientes. Los IDs y los identificadores de color se mostrarán en el diseño de la placa.
Eliminar selección	Solo en el modo Diseño de placas. Este comando borrará los IDs y los identificadores de colores de los pocillos seleccionados y los dejará en blanco.
Seleccionar todas las placas sin usar	Solo en el modo Diseño de placas. Se marcan todos los pocillos sin usar de la placa.
Crear o eliminar alias	Solo en el modo Diseño de placas. Establece o elimina designaciones de alias para los nombres de los pocillos. Consulte la sección 4.3.2 Diseño de método: Cómo definir un diseño de placas/ Asignar alias al pocillo requerido.

Barra de control de la pestaña Crear o editar método

La **Barra de control** a la izquierda de la pantalla ofrece un número de opciones que deberían ejecutarse en la secuencia sugerida. En función del tipo de medición y el tipo de lector conectado, algunas opciones puede que no estén disponibles y estén ocultas.

Al seleccionar un elemento en la barra de control se mostrará el cuadro o la vista de placa correspondiente en el panel de **Ventana de diseño de placas** a la derecha de la pantalla.

Todas las opciones disponibles se marcan con una marca de control, una vez se han definido. Para cualquier tipo de transformación aparece el nombre de transformación.

La pestaña **Barra de control - Crear o editar método** contiene los elementos siguientes:



Grupo	Elemento Diseño de placas
Diseño de método	La ventana Diseño de placas se abre, mostrando el diseño de placas y el cuadro de diálogo Asignación de los pocillos .
	Elemento Valores Conc., Dil. y Ref.
	Se abre el cuadro de diálogo de definición de Concentración, dilución o referencia y muestra el valor de dilución, concentración o referencia de cada pocillo. Se pueden establecer las concentraciones estándar, así como los factores de dilución y los valores de referencia. Una función de Llenado automático ofrece una asignación fácil de las concentraciones en caso de relaciones matemáticas peculiares de las concentraciones entre los pocillos individuales.
Cálculo previo	 Reducción de datos de espectros para exploraciones de longitud de onda
Grupo Datos transformados	 Elemento Añadir nueva transformación Se abre, la ventana Diseño de placas y muestra el diseño de la placa y el campo de entrada del diálogo Transformaciones.
Grupo Cinético	 Elemento Reducción de datos cinéticos Se abre la ventana Parámetros de cálculos cinéticos. Se puede definir la evaluación de los datos de medición cinética.
Transformación cinética	 Elemento Añadir nueva transformación cinética Se abre la ventana Diseño de placas y muestra el diseño de la placa y el campo de entrada del diálogo Transformaciones (solo se pueden seleccionar datos de entrada cinéticos).
Grupo Concentraciones	 Elemento Curva estándar La ventana Curva estándar se abre para fijar los parámetros para el cálculo de concentraciones y la presentación gráfica de la curva estándar de los datos evaluados.
Grupo Datos de transformación de concentración	 Elemento Añadir nueva transformación de concentración Se abre la ventana Diseño de placas y muestra el diseño de placa y el campo de entrada del diálogo Transformaciones de concentración (como datos de entrada, solo se pueden seleccionar las concentraciones).
Grupo Evaluar datos	 Elemento Definición del corte Elemento Validación QC Se abre la ventana Definir cortes. En este cuadro de diálogo se pueden definir los límites para una evaluación cualitativa (selección). Se abre la ventana Definir validación QC. Se utiliza para verificar la validez de una prueba.
Grupo Procesamiento de datos	 Elemento Exportación de datos Elemento Informe impreso Elemento Procesamiento automatizado de datos



Grupo Otras opciones	Elemento Mensajes al usuario
	Elemento Formato de números
	Elemento Notas de método
	La ventana Definir mensajes al usuario permite asignar datos (palabras clave, comentarios o mensajes) a cada medición, que luego pueden incorporarse en una impresión.
	La ventana Formato de números permite al usuario definir el formato de los números para los datos sin procesar o los datos transformados que se muestran.
	En Notas de método se puede escribir una descripción del método.

Barra de control - Pestaña Crear o editar método: menú contextual

Haciendo clic con el botón secundario del ratón en cualquier transformación, transformación cinética y transformación de concentración en la barra de control, aparece un **menú contextual** para la transformación marcada. Están disponibles los siguientes comandos:

Menú	Descripción
Cambiar nombre de transformación	Se puede asignar un nombre distinto a la transformación seleccionada.
Insertar transformación	Se utiliza para definir una nueva transformación.
Quitar transformación	Se utiliza para borrar una transformación.



4.3.2 Diseño de método: Cómo definir un diseño de placas

En la Barra de control expanda Diseño de método y seleccione Diseño de placas. Aparecen la Vista de placa y el cuadro de diálogo Asignación de los pocillos.

Para cada pocillo en la placa de análisis se puede asignar un identificador. En base a esto debe definirse el diseño de placas, (p. ej. a un pocillo que deba aparecer como un control positivo debe asignársele un identificador diferente que a otro pocillo que deba aparecer como un control negativo).

Los identificadores estándar son:

Asignación de los pocillos	Muestra	SM (muestra)
Identificador Grupo exp.:	Blanco	BL (blanco) BF (búfer de referencia de polarización)
Núm. de ID.	Referencia	RF (referencia de polarización)
	Estándar	ST (estándar)
NC LPC HPC BF RF Def. identif.	Control	PC (control positivo) NC (control negativo) LPC (control positivo bajo) HPC (control positivo alto) CL (calibrador)
Selección de Borrar relleno selección		

Asignar un identificador al pocillo requerido

En el cuadro de diálogo **Asignación de los pocillos** se puede realizar la identificación de los pocillos y la definición del diseño. Este cuadro también ofrece un número de capacidades de asignación de ID automatizadas, que son una herramienta esencial para placas de alta densidad.

La configuración deseada debe realizarse en el cuadro de diálogo **Asignación de los pocillos**. Generalmente, los pocillos se pueden seleccionar haciendo clic en el pocillo individual o arrastrando el ratón sobre los pocillos necesarios.

Existen las siguientes maneras de asignar el identificador seleccionado a los pocillos:

- Haga doble clic al realizar la selección del pocillo
- Seleccione los pocillos de la microplaca y luego haga clic en el botón Rellenar selección (o haga clic con el botón secundario del ratón y seleccione Rellenar selección en el menú contextual) en el cuadro de diálogo Asignación de los pocillos
- Seleccione los pocillos en la microplaca y luego haga doble clic en el identificador en el cuadro de lista de identificadores del diálogo de asignación de los pocillos.

Tras asignar las definiciones cambiará la pantalla en los pocillos modificados.

SM1_4	1ª línea: muestra, número de grupo experimental de 1, número de ID de muestra de 4.	
1/14	2ª línea: número de réplicas es 1, número total de réplicas es 14.	
x-BL1	3ª línea: vacía en principio – rellenada con la fórmula de transformación o un valor de concentración, dilución o referencia si se define, p. ej. reducir el valor de este pocillo (x) mediante los blancos.	
El cuadro de diálogo Asignación de los pocillos contiene los siguientes elementos:		
Cuadro de grupo	Seleccione identificadores de la lista desplegable para	

Ejemplo de un pocillo etiquetado:

elementos:	
Cuadro de grupo Identificadores	Seleccione identificadores de la lista desplegable para marcar los pocillos.
	 Haga clic en el botón Definir identif para definir identificadores adicionales.
Campo de selección Grupo exp.	Si la placa está formada por más de una prueba, p. ej. más de un grupo experimental, defina el Grupo experimental al que pertenecen los pocillos.
Cuadro de grupo Núm. de ID	 El Número de ID se utiliza para asignar el mismo ID a las réplicas que van juntas. El número de ID solo está disponible para muestras y estándares. Campo de selección Núm. de ID La opción Núm. de ID permite al usuario marcar aquellas réplicas que se asocian con el mismo ID. Esto se puede usar si las réplicas se han asignado a diferentes áreas de la placa. Botones de flecha El ID se calcula automáticamente. Si se marca un grupo de pocillos, mediante los botones de flecha se puede establecer la dirección en que se asignan los IDs a los pocillos (verticalmente, horizontalmente).
Cuadro de grupo Réplicas	 Determina el número de réplicas para el tipo de identificador seleccionado: Botón de opción Número fijo Solo activado para estándares y muestras donde se puedan usar IDs. Este número define cuántas réplicas se destinan a este identificador. Botón de opción Todo Todos los pocillos seleccionados se definen como réplicas de un identificador. Si se elige un número de ID existente para las muestras y estándares, los pocillos seleccionados se añaden como réplicas de las réplicas existentes. Con el resto de tipos de identificadores los pocillos se añaden como réplicas de las réplicas existentes. Dos Botones de flechas definen la dirección para contar el número de réplica.
Botón Rellenar sección	Si se ha marcado un área de la placa, puede rellenarse con los identificadores correspondientes. Los IDs y el color de los identificadores se mostrarán en el diseño de las placas.



Botón Eliminar sección	Haga clic en Eliminar o pulse SUPR para borrar los IDs y los identificadores de colores de los pocillos seleccionados y dejarlos en blanco.
Botón	Haga clic en Def. identif . si hay que definir un identificador
Definir	nuevo o si hay que editar un identificador existente (consulte
identificador	Definir un nuevo identificador)

Asignar alias al pocillo requerido

Para asignar denominaciones de alias para nombres de pocillos definidos, haga clic con el botón secundario del ratón en el pocillo requerido, seleccione **Crear o quitar alias...** y seleccione un nombre de identificador alternativo en la lista desplegable. El alias se marca con un asterisco * y tiene el mismo grupo experimental, número de ID y número de réplica que el pocillo definido primariamente. Esta función se utiliza, p. ej. si el estándar 0 también se utiliza como control negativo.

Definir un nuevo identificador

En el cuadro de diálogo **Asignación de pocillos** haga clic en **Def. identif.** para definir un nuevo identificador o editar uno existente.



Los identificadores se enumeran en grupos (véase la tabla siguiente). Al resaltar un identificador, sus propiedades se muestran en la ventana de la derecha.

Los identificadores estándar son:

Muestra	SM (muestra)
Blanco	BL (blanco)
	BF (búfer de referencia de polarización)
Referencia	RF (referencia de polarización)
Estándar	ST (estándar)
Control	PC (control positivo)
	NC (control negativo)
	LPC (control positivo bajo)
	HPC (control positivo alto)
	CL (calibrador)

El cuadro de diálogo Definir identificadores contiene los siguientes elementos:



Estructura de árbol Identificador	En una pequeña ventana se ofrece una vista estructurada de todos los identificadores existentes, sus colores y abreviaturas. Los identificadores se enumeran bajo los grupos Muestra, Blanco, Referencia, Estándar y Control .
Cuadro de grupo Identificador	Se mostrarán los criterios mantenidos por los diversos identificadores utilizados en el programa. Si se necesitan nuevos identificadores, pueden insertarse aquí. En el cuadro de grupo Identificador se muestra el tipo, la abreviatura, el color del pocillo y la descripción del identificador seleccionado en la estructura de árbol
	El cuadro de grupo Identificador solo se activa para introducir datos cuando
	 se selecciona uno de los grupos de identificadores en la estructura de árbol y se hace clic en el botón Insertar, o bien
	 se selecciona uno de los Identificadores en la estructura de árbol y se hace clic en el botón Editar.
	Entonces, en este cuadro de grupo será preciso definir el tipo, la abreviatura, el color del pocillo y la descripción del nuevo identificador.
	Campo de texto Tipo : Muestra el tipo del identificador. No permite cambios
	Campo de texto Abreviatura : Se usará la abreviatura del identificador en la pantalla de placa de análisis.
	Lista desplegable Color : Aquí se debe seleccionar el color del identificador para el diseño de placa.
	Campo de texto Descripción : Para cada identificador se puede introducir una descripción de texto.
	Use uno de estos dos botones para guardar los cambios de datos recién introducidos o desechar los cambios:
	Botón Guardar : El botón Guardar guarda el color, la abreviatura y la descripción que se han introducido para el identificador correspondiente.
	Botón Desechar: El botón Desechar cancelará cualquier cambio.
Botón Insertar	Haga clic en el botón Insertar para crear un nuevo identificador. Este nuevo identificador se asociará con el grupo de identificadores seleccionado actualmente en la estructura de árbol.
Botón Editar	Los identificadores seleccionados se pueden modificar.
Botón Eliminar	Haga clic en Eliminar o pulse Supr para borrar un identificador seleccionado.



Botón Establecer como predeterminado	Usando esta opción se pueden definir los ajustes como predeterminados para su uso futuro.
Botón Restablecer valores predeterminados	Usando esta opción se pueden restablecer los ajustes a los valores predeterminados definidos previamente.

4.3.3 Diseño de método: Valores Conc., Dil. y Ref.

En la Barra de control expanda Diseño de método y seleccione Valores conc., dil. o ref. Aparecen la Vista de placa y el cuadro de diálogo Seleccionar identificador.



La ventana **Concentración, dilución o referencia** contiene los elementos siguientes:

Lista desplegable Seleccionar identificador	Los valores están correlacionados con los tipos de pocillos individuales; este campo ofrece todos los pocillos actualmente incluidos en la medición. Seleccione un identificador para editar la dilución correspondiente o los factores de concentración. Se listan todos los pocillos que coinciden con el identificador del tipo de pocillo seleccionado.
Campo de selección Grupo exp.	Seleccione el grupo experimental correspondiente. Si solo existe un grupo experimental, el campo no se puede editar.



Tabla con columnas Identificador y Concentración, dilución	Esta tabla muestra todos los pocillos en la columna Identificador que coinciden con el identificador seleccionado (a partir de la lista Seleccionar identificador). Los valores de dilución, concentración o referencia correspondientes se insertan y se enumeran en la columna Dilución, concentración o referencia. P. ej. introduciendo un factor de dilución de 2 significa que la muestra se ha diluido a la mitad. Por tanto, la concentración calculada se multiplica por 2.
Campo de texto Unidad	Se puede determinar la unidad de concentración mostrada.
Casilla de verificación Llenado automático	La función Llenado automático ofrece el cálculo de las concentraciones correspondientes o factores de dilución de acuerdo con los tipos de series disponibles.
Botones Opción para el cálculo matemático de concentración	 Si se ha seleccionado la casilla de verificación Llenado automático, existen las opciones siguientes: Series aritméticas: a (n+1) = a (n) + Series geométricas: a (n+1) = a (n) x Series definidas por el usuario a(n+1) = Ejemplo: Cada concentración posterior debería tener el doble + 0,5 de la concentración anterior: Fórmula 1: a(n+1) = 2 * n + 0,5
Botón Aplicar	El botón Aplicar aplica el cálculo matemático de concentración seleccionado a los pocillos como se muestra en la tabla con columnas Identificador y Concentración, dilución .
Cuadro de grupo Series de dilución	Casilla de verificación Calcular ICx Muestra un gráfico de dilución de la muestra y calcula automáticamente los valores ICx. Esto requiere muestras con al menos cuatro réplicas y al menos cuatro diluciones diferentes definidas*.
	Lista desplegable Datos de entrada Seleccione de la lista desplegable los datos de entrada.
	Campo de texto Condición de cálculo La intercepción se calcula con el porcentaje introducido del valor máximo en relación al valor máximo menos el valor mínimo*.
	Campo de texto Nombre de ICx Se pueden seleccionar datos para el cálculo ICx. El nombre se rellena automáticamente en función de la condición de cálculo.
	Establecer valor de 0% como Intensidad 0* Intensidad mín. de series de dilución*

La dilución se puede editar directamente en la ventana de diseño de placas haciendo clic con el botón secundario en uno o más pocillos seleccionados. De este modo es posible asignar diferentes valores de dilución a las réplicas.



Cálculo ICx

El cálculo matemático del ajuste para las series de dilución es idéntico al cálculo de la curva estándar con el algoritmo de Marquardt con 4 parámetros.

Requiere al menos cuatro réplicas con distintas diluciones.

Además, se calculan las intercepciones especificadas, p. ej., IC 50.

Destaced Reference Property	10% 100% Zoom														- î
Decis de mitodo Decis de mitodo Decis de las placas Valoras de Conc., DL., Raf	Selectors: 24 Graditator: 1 1 1 1 1 2 2 2 2 3 3 3		1	2	3	4,	5	6	7	8	9	10	11	12	
Allade nueves transformació	4 4 4 4 4 4 6 6 7 7 7 7 7 7 7 7 7 7 7 7 7 7 7 9 9 9 9 9 9 9 9 9 9 9 9 9 9 9 10 10 10 10 10 10 10 10 10 10 10 10 10 10 10 10 10 10 10 10 10 10 10 10 10 10 10 10 10 10 10 10 10 10 10 10 10 10 10 10 10 10 10 10 10 10 10 10 10 10 10 10 10 10 10 10 10 10 10 10 10 10 10 10 10 1	A	ST1_1 1/1 0	SM1_1 1/11 1	SM1_1 2/11 2	SM1_1 3/11 3	SM1_1 4/11 4	SM1_1 5/11 5	SM1_1 6/11 6				SM1_1 10/11 10	SM1_1 11/11 11	
Afade nue la transformació Note transformació Definición del carte Validación QC	11 H		ST1_2 1/1 1	SM1_2 1/11 1	SM1_2 2/11 2	SM1_2 3/11 3	SM1_2 4/11 4	SM1_2 5/11 5	SM1_2 6/11 6	SM1_2 7/11 7	SM1_2 8/11 8	SM1_2 9/11 9	SM1_2 10/11 10	SM1_2 11/11 11	
Procesaniesto de d		C	ST1_3 1/1 2	SM1_3 1/11 1	SM1_3 2/11 2	SM1_3 3/11 3	SM1_3 4/11 4	SM1_3 5/11 5	SM1_3 6/11 6	SM1_3 7/11 7	SM1_3 8/11 8	SM1_3 9/11 9	SM1_3 10/11 10	SM1_3 11/11 11	
Otras opciones S Indicaciones al casario Pornato de números Notas de reétado			ST1_4 1/1 3	SM1_4 1/11 1	SM1_4 2/11 2	SM1_4 3/11 3	SM1_4 4/11 4	SM1_4 5/11 5	SM1_4 6/11 6	SM1_4 7/11 7	SM1_4 8/11 8	SM1_4 9/11 9	SM1_4 10/11 10	SM1_4 11/11 11	
	vedati ngt			SM1_5 1/11 1	SM1_5 2/11 2	SM1_5 3/11 3	SM1_5 4/11 4	SM1_5 5/11 5	SM1_5 6/11 6	SM1_5 7/11 7	SM1_5 8/11 8	SM1_5 9/11 9	SM1_5 10/11 10	SM1_5 11/11 11	
	- a(n+1) - a(n) + 0 = a(n+1) - a(n) × 0 = a(n+1) -	F	ST1_6 1/1 5	SM1_6 1/11 1	SM1_6 2/11 2	SM1_6 3/11 3	SM1_6 4/11 4	SM1_6 5/11 5	SM1_6 6/11 6	SM1_6 7/11 7	SM1_6 8/11 8	SM1_6 9/11 9	SM1_6 10/11 10	SM1_6 11/11 11	m
	Series de diución Catolier ICx Dotos en proceser Catolic cond. (%) Rombre			SM1_7 1/11 1	SM1_7 2/11 2	SM1_7 3/11 3	SM1_7 4/11 4	SM1_7 5/11 5	SM1_7 6/11 6	SM1_7 7/11 7	SM1_7 8/11 8	SM1_7 9/11 9	SM1_7 10/11 10	SM1_7 11/11 11	CAN Austri
	1 2 2 2 2 2 2 2 2 2 2 2 2 2 2 2 2 2 2 2	Н	ST1_8 1/1 7	SM1_8 1/11 1	SM1_8 2/11 2	SM1_8 3/11 3	SM1_8 4/11 4	SM1_8 5/11 5	SM1_8 6/11 6	SM1_8 7/11 7	SM1_8 8/11 8	SM1_8 9/11 9	SM1_8 10/11 10	SM1_8 11/11 11	TE
	O promocio min. de serie de dikudin														

Es posible definir más de una intercepción para las series de dilución.

Para definir el valor de 0% para el cálculo, seleccionar:

- Intensidad 0
- Usar 0 DO

El mayor valor de las series de dilución se considera el 100%, el valor de 0 se considera 0%. El ICx (p. ej., IC 50) entonces se define como la dilución en la que la respuesta alcanza el x% (p. ej., el 50%). Solamente se calcula si el valor está dentro de los datos disponibles (no se utiliza ninguna extrapolación).

O bien,

Intensidad mínima de las series de dilución
 Utilizando la intensidad mínima de las series de dilución



Aviso

El IC 50 a menudo se calcula con el valor medio de dilución=1:1 y dilución=1:infinito. Para conseguirlo, las series de dilución se deben corregir según su línea base antes de calcular ICx.

El mayor valor de las series de dilución se considera el 100%, la intensidad mínima de las series de dilución se considera el 0%.



4.3.4 Cálculo previo: Reducción de datos de espectros

Esta opción solo está disponible para mediciones que contengan una exploración bidimensional (exploración de absorbencia).

Expanda Cálculo previo en la barra de control y seleccione Reducción de datos de espectros.

La reducción de datos de espectros permite realizar una reducción del espacio en blanco de espectros y extraer datos específicos, p. ej. intensidades, de la exploración.

La ventana **Reducción de datos de espectros** contiene los elementos siguientes:

Reducción del blanco	Opciones	Nembro do datos
Suavizar	Opciones	S Espectros suavizados
Selección de longitud de onda	Opciones	S Espectro normalizado
Normalizar	Opciones	V Pico de longitud de onda
✓ Pico	Opciones	V Pico de anchura V Pico de área
Fórmulas personalizadas	Opciones	

Casilla de verificación Reducción del blanco	Esta opción solamente está disponible si hay blancos definidos en el diseño. Los espectros reducidos blancos se calculan restando el espectro del pocillo de blanco de todos los demás pocillos. Cuando hay más blancos definidos en la placa, p. ej., cuando se definen varios grupos experimentales, se muestra un botón de opciones adicional que permite definir qué blanco se debe utilizar.
	Cuadro de texto Datos de entrada : muestra los datos de entrada que se procesarán.
	Botón de opción Reducción del blanco : seleccione si desea efectuar una reducción del blanco en el grupo experimental seleccionado.
	Lista desplegable Reducir todos por : seleccione para definir el identificador que debe utilizarse para la reducción del blanco.
Casilla de verificación Suavizar	Permite suavizar espectros que contienen mucho ruido. En la sección de opciones se puede configurar un factor de suavizado. Ese factor define el grado de suavizado.
	Cuadro de texto Datos de entrada : muestra los datos de entrada que se procesarán.
	Factor de suavidad: definir el factor de suavizado.



Casilla de verificación Selección de longitud de onda **Selección de longitud de onda** se usa para extraer intensidades a longitudes de onda específicas y calcular resultados como relaciones. Es obligatorio establecer las opciones.

El cuadro de diálogo de parámetros de **Selección de longitudes de onda** contiene los siguientes elementos:

Parámetros d	e selección de lo	ongitudes de o	nda			×
Entrada:	Espectros suav	izados				
 Selecciona Calcular p Área A definir p 	ar longitudes de o roporción por el usuario Longit	nda WL [nm] WL [nm] de WL [nm] Fórmula udes de onda 500	500 600 500 Constantes:	/ WL [nr para WL [nr] 700 i] 700 Dperadores + v	>
Nombre de 500 nm 600 nm / 70 <	datos 00 nm Cancelar	Formula x[500] x[600]/x[700	0]		>	Anadir Borrar Editar Ayuda
Se puede específic pueden i	en extrae a, se pue ntroducir	r intensic den calc fórmulas	lades a ular pro person	una longi porciones alizadas.	tud de o s y área	onda s, y se
Cuadro o entrada o	de texto D que se pro	atos de	entrada	a: muestra	a los da	tos de
Botón de puede af intensida	e opción S ňadir a la id a una lo	Seleccio lista de r ongitud c	nar long esultado le onda	jitudes d os calcula específic	e onda ados una a.	: se a
Botón de especific entre inte muestra	e opción C ar dos lor ensidades el resulta	alcular ngitudes a esas do.	propore de onda ongitud	ción : perr a. Se calc es de onc	nite ula la re da y se	elación
Área: pa entre dos	ra calcula s longitud	ar el área es de on	bajo la da defin	curva de iidas.	especti	OS
Botón de introduci intensida pueden u	e opción A r una fórm ides a lon usar en fó	definir nula defir gitudes o rmulas.	por el ι nida por de onda	i suario : p el usuari especific	bermite o. Las adas se)
Cuadro o definidas	de lista Re s.	esultado	s : resur	ne todas	las fórm	nulas
Botón Aí resultado	ñ adir : la s os.	elección	actual	se añade	a la list	a de
Botón Bo	orrar: el re	esultado	seleccio	onado se	borra la	a lista.
Botón Ec del resul	litar : se p tado selec	ouede de ccionado	finir el n	ombre pa	ara los c	latos



Casilla de verificación	El cuadro de diálogo Normalizar parámetros contiene los siguientes elementos:								
Normalizar	Normalizar parámetros X								
	Entrada: Espectros suavizados OK								
	Normalizar a Cancelar								
	Intensidad max								
	O Intensidad con LO [nm] 400 Ayuda								
	Utilice esta opción para normalizar el espectro a la intensidad a una longitud de onda dada o a la intensidad máxima del gráfico de espectro de cada pocillo.								
	Cuadro de texto Datos de entrada : muestra los datos de entrada que se procesarán.								
	Normalizar a: seleccione entre intensidad máxima o inserte una intensidad personalizada.								
Casilla de verificación Pico	Un pico se define como la posición del valor máximo del espectro. Los desbordamientos y los máximos en el borde o izquierdo o derecho del espectro no se tratan como picos. Es muy importante suavizar suficientemente el espectro antes de buscar un pico.								
	siguientes elementos:								
	Pico de parámetros X								
	Entrada: Espectro normalizado OK								
	Rango de De: 400 nm Cancelar								
	A: 600 nm								
	Omitir picos interiores a 0 OD/RFU								
	Calcular anchura y área en 50 % de intensidad máx Ayuda								
	Utilice esta opción para encontrar el pico con la intensidad más alta dentro de una longitud de onda definida.								
	También se pueden definir el umbral (valor DO/RFU) y los criterios de cálculo para calcular la anchura y el área.								
	Se calcularán los datos siguientes para la ruta encontrada:								
	pico de intensidad								

- longitud de onda del pico
- anchura del pico



Casilla de verificación Fórmulas	Seleccione esta casilla para introducir fórmulas para calcular el espectro utilizando las funciones dadas; p. ej. suavizar.					
personalizadas	Para obtener información adicional, consulte la sección 11.2 Reducción de datos de espectros.					
	La transformación debe devolver un espectro o un dato único para todos los pocillos; no es posible mezclar ambos. Utilizando esta opción se pueden definir funciones que no están disponibles en la selección estándar. P. ej., es posible encontrar más picos dentro de un espectro o crear una variación de un espectro.					
Resultados disponibles	Se enumeran los resultados disponibles de todos los datos. S es un espectro					
	v es un valor.					

4.3.5 Datos transformados: Añadir nueva transformación

En la barra de control expanda **Datos transformados**. Todas las transformaciones definidas se muestran en la barra de control.

Para definir una nueva transformación, seleccione los pocillos a los que se debe aplicar la transformación y haga clic en **Añadir nueva transformación...**.

El nombre predeterminado debe cambiarse inmediatamente en el cuadro de edición. El nombre también puede insertarse o cambiarse más tarde haciendo clic en **Cambiar nombre de transformación** desde el menú contextual.

Aviso

Los nombres de las transformaciones deben usar exclusivamente caracteres latinos para que dichas transformaciones estén disponibles como datos de entrada adicionales para transformaciones posteriores con datos de entrada múltiples.



Aviso

El nombre de una transformación se utiliza para representar el resultado de los cálculos y también se utiliza cuando se muestran los valores en los menús de emisión de datos específicos. Los valores calculados de las transformaciones también están disponibles como datos de entrada para otras evaluaciones.

Un ejemplo típico de una transformación es la reducción del valor vacío (espacio en blanco) de todos los pocillos.



Si se fija un espacio en blanco en el Diseño de placas, la fórmula para calcular la reducción del espacio en blanco se ofrece en el cuadro combinado: x-BL1.

El símbolo x hace referencia al valor actual en un pocillo. BL1 es el valor medio de los pocillos en blanco.





Aviso

Si se define una medición de etiquetas múltiples con dos etiquetas, en el cuadro combinado se ofrecen diferentes fórmulas para reducir los datos de etiquetas múltiples.

'Etiqueta1'!x/'Etiqueta2'!x ... cálculo de relación predefinida 'Etiqueta2'!x/'Etiqueta1'!x ... cálculo de relación predefinida 'Etiqueta1'!x-'Etiqueta2'!x ... cálculo de diferencia predefinida 'Etiqueta2'!x-'Etiqueta1'!x ... cálculo de diferencia predefinida ('Etiqueta2'!x-'Etiqueta1'!BL1)/('Etiqueta2'!x-'Etiqueta2'!BL1) ... cálculo de relación predefinida con reducción del blanco ('Etiqueta2'!x-'Etiqueta2'!BL1)/('Etiqueta1'!x-'Etiqueta1'!BL1) ... cálculo de relación predefinida con reducción del blanco



Aviso

Si el modo de lectura se ha establecido como absorbencia, en el cuadro combinado se mostrará la fórmula para calcular la transmisión: 1/10^x.

Al configurar el método se pueden definir varios cálculos de transformación. Un ejemplo típico de ello es la deducción del valor vacío (blanco) de todos los pocillos. Para obtener detalles adicionales, consulte la sección 11.3 Escribir una fórmula.

El cuadro de texto Fórmula encima de la vista de placa se usa para definir estos varios cálculos de transformación utilizando datos sin procesar y transformaciones anteriores como datos de entrada.

En la microplaca se deben seleccionar los pocillos a los cuales se va a aplicar la transformación.

El campo de edición de transformaciones contiene los siguientes elementos:

Lista desplegable Entrada	Esta lista contiene los valores de mediciones, los resultados de los cálculos previos y todas las transformaciones ya definidas, así como los promedios. Si se seleccionan datos Media como datos de Entrada para Transformaciones , entonces se asignarán transformaciones a la primera réplica de un identificador solamente
Cuadro de texto Fórmula fx	Las fórmulas se pueden introducir escribiéndolas o seleccionando la función requerida en el cuadro de lista Funciones&Constantes.
Lista desplegable Fórmula	Esta lista contiene algunas fórmulas estándar (véase las notas a continuación) y todas las fórmulas que se han incorporado al método actual. Por lo tanto, se puede seleccionar una fórmula en esta lista o se puede añadir una nueva fórmula. Se pueden introducir fórmulas para transformaciones utilizando los correspondientes operadores, variables y funciones.
Botón Confirmar	Modo de edición de fórmula solamente. Marca de verificación verde. Asigna las transformaciones al pocillo seleccionado. Haga clic en el botón Confirmar o pulse la tecla INTRO para asignar la definición de fórmula al pocillo y cambiar al modo Seleccionar .



Botón Cancelar	Modo de edición de fórmula solamente. Aspa roja. Haga clic en el botón Cancelar o pulse la tecla CANCELAR para salir del modo Editar sin asignar la definición de fórmula al pocillo.
Lista desplegable Datos disponibles	Seleccione en esta lista los datos apropiados si piensa usar para los cálculos más de un conjunto de datos de entrada. El conjunto de datos se mostrará entre apóstrofos seguido de un signo de admiración en el cuadro de texto Fórmula. Para completar la definición, introduzca el nombre del identificador o consulte el valor correspondiente dentro del pocillo. P. ej., 'Datos sin procesar' ! BL1
	La lista contiene los valores de mediciones, los resultados de los cálculos previos y todas las transformaciones definidas hasta ese momento.
Lista desplegable Funciones	Enumera todas las funciones matemáticas y booleanas para la definición de fórmulas.
Botón Opciones	 Utilice este botón para personalizar el comportamiento cuando seleccione pocillos en el modo de edición. Puede seleccionar entre las siguientes opciones: Nombre de media de identificador Nombre de réplica de identificador Nombre de pocillo de placa
Botón Constantes	Haga clic en este botón para abrir el cuadro de diálogo Definir constantes para definir constantes para los cálculos.

La definición de transformación tiene dos modos:

Modo Seleccionar	Al hacer clic en un pocillo, en el cuadro de texto se muestra la fórmula correspondiente. El pocillo activo en ese momento se muestra con un borde rojo.
Modo Editar	Cuando se introduce una fórmula o se pulsa la tecla "=", el pocillo activo se muestra con un borde azul. Al seleccionar otros pocillos se agrega el identificador correspondiente a la fórmula actual. Cuando se selecciona el pocillo activo, se añade una "x" a la fórmula. El símbolo "x" se refiere al valor actual dentro de un pocillo.

Una vez que se crea la fórmula, se asignará al pocillo seleccionado al pulsar Intro o al hacer clic en el botón de confirmar (marca de verificación verde). Después de asignar la fórmula, se accede al modo seleccionar. Un comprobador de fórmulas integrado verifica la plausibilidad de la fórmula. En caso de que la fórmula creada no sea ejecutable, se muestra un aviso.

Seleccione **CANCELAR** para desechar la fórmula definida y acceder al modo seleccionar. También se pueden utilizar los botones correspondientes a la izquierda del cuadro de texto Fórmula.

Cuando esté en modo **Seleccionar**, haga clic en el cuadrado rojo del borde inferior derecho del pocillo y arrastre el marco de selección rojo sobre los pocillos a los que desea aplicar la fórmula.

También puede seleccionar los pocillos deseados y utilizar las funciones **Crear fórmula** y **Quitar fórmula(s)** disponibles en el menú contextual.



Estas acciones se deben realizar para todos los pocillos, que contendrán transformaciones.

Definir constantes

Utilice el cuadro de diálogo **Definir constantes** para definir los valores constantes de un método. Estas constantes se pueden utilizar dondequiera que se pueda escribir una fórmula.

Haga clic en el botón Aceptar para guardar los parámetros editados.

El cuadro de diálogo Definir constantes contiene los siguientes elementos:

Lista	En cada fila se puede definir una constante. Las filas están
Constantes	divididas en cuatro columnas:

Nombre

Escriba un nombre de identificador apropiado para la constante, por ejemplo: un código o una abreviatura. Solamente se pueden utilizar letras.

Valor

En el campo de texto **Valor** se debe asignar un valor numérico a la constante.

Comentario

Escriba un comentario corto relativo a la constante.

Req. ("requerida")
 La casilla de verificación Req. indica que una medición

solamente se puede ejecutar al recibir o confirmar el valor para la constante.

4.3.6 Cinética: Reducción de datos cinéticos

En la barra de control expanda **Cinética** y seleccione **Reducción de datos cinéticos**.

El cuadro de diálogo está estructurado en pestañas: **Pendientes**, **Iniciaciones**, **Mín./Máx./Área**, **Datos de salida disponibles**, **Cinética de enzima**. Las unidades de los valores se muestran según el modo de medición seleccionado (por ejemplo: OD para absorbencia).



Aviso En el cálculo de datos cinéticos, no se tienen en cuenta los datos inservibles (p. ej., los valores de desbordamiento).

Pestaña Pendientes

Esta pestaña permite al usuario establecer la evaluación de la pendiente de la curva cinética:

Lista desplegable Datos de entrada	Seleccione los datos de entrada que desea procesar.
Lista desplegable Cálculo	Seleccione el método de cálculo, lineal o cuadrático (consulte el capítulo 11 Cálculos).
Casilla de verificación Pendiente media	En los campos de texto suministrados se pueden indicar los tiempos de inicio y fin, a no ser que se desee analizar toda la cinética.



	 Botones de opción Tiempo/Puntos: si se selecciona Tiempo, se deben escribir los tiempos de inicio y fin en horas, minutos y segundos. si se selecciona Puntos, se debe escribir el número de ciclos cinéticos para indicar los puntos de inicio y de final del análisis.
	 Campo Inicio: aquí se debe escribir la hora de comienzo (en horas, minutos y segundos) o el número de ciclos.
	 Campo Final: aquí se debe escribir la hora de finalización (en horas, minutos y segundos) o el número de ciclos.
Casilla de verificación Pendiente máxima	En los campos de texto suministrados se pueden indicar los tiempos de inicio y fin, a no ser que se desee analizar toda la cinética.
	 Botones de opción Tiempo/Puntos: si se selecciona Tiempo, se deben escribir los tiempos de inicio y fin en horas, minutos y segundos. si se selecciona Puntos, se debe escribir el número de ciclos cinéticos para indicar los puntos de inicio y de final del análisis.
	 Campo Inicio: aquí se debe escribir la hora de comienzo (en horas, minutos y segundos) o el número de ciclos.
	 Campo Final: aquí se debe escribir la hora de finalización (en horas, minutos y segundos) o el número de ciclos.
	 Campo de texto Puntos: seleccione para cuántos puntos se realiza el cálculo de la pendiente máxima.

La opción **pendiente media** determina la pendiente media (aumento/disminución media) a lo largo del intervalo definido. Primero se deben introducir los tiempos de inicio y de finalización y un modo de cálculo (lineal o cuadrático). Se determinarán todos los puntos de medición que están comprendidos dentro del intervalo seleccionado. Se trazará una línea de regresión (regresión lineal o polinomio de 2º grado) a través de los puntos de medición seleccionados y se conformará la pendiente media. La pendiente media se define como la media aritmética de las pendientes que se calculan desde los puntos centrales de dos puntos de medición adyacentes. Los datos de cinéticas accesibles mediante este método incluyen la pendiente media por segundo, por minuto y por hora, así como el coeficiente de correlación y la calidad del ajuste.

La pendiente máxima (aumento/disminución máxima) conseguida en el intervalo seleccionado se determinará con la opción **pendiente máxima**. En la opción **Puntos** es preciso introducir el número de puntos combinados. Para empezar, la pendiente se calculará desde el punto central del 1.º y el n.º punto desde los primeros n puntos. Después el intervalo se moverá un punto adelante y se repetirá el proceso. Ese proceso continuará sobre todos los puntos dentro del intervalo seleccionado. El resultado se determinará como el mayor valor absoluto de estas pendientes individuales. Los datos de cinéticas accesibles mediante este método incluyen la pendiente máxima por segundo, por minuto y por hora, así como el lapso, en segundos, desde la primera medición hasta la pendiente máxima.



Pestaña Iniciaciones

Utilice la pestaña **Iniciaciones** para determinar el tiempo hasta alcanzar un cierto punto de datos (valor de DO):

Lista desplegable Datos de entrada	Seleccione los datos de entrada que desea procesar.
Casilla de verificación Tiempo hasta iniciación	Si está activada la casilla de verificación Tiempo hasta iniciación , se puede escribir en el siguiente campo de texto un valor absoluto para el comienzo.
	 Campo de texto: se debe escribir un valor absoluto para la iniciación.
	 Botón de opción y campo de texto Base media de los primeros n puntos: si lo selecciona, escriba el número de puntos que desee.
	 Botón de opción y campo de texto Base: si lo selecciona, debe escribir en el campo de texto adyacente un valor absoluto para la base.
Casilla de verificación Tiempo hasta iniciación %	Si está activada la casilla de verificación Tiempo hasta iniciación %, se puede escribir en el siguiente campo de texto un valor de porcentaje.
	 Campo de texto: se debe escribir un valor de porcentaje para la iniciación.
	 Botón de opción y campo de texto Base media de los primeros n puntos: Si lo selecciona, puede escribir el número de puntos que desee.
	 Botón de opción y campo de texto Base: si lo selecciona, debe escribir en el campo de texto adyacente un valor absoluto para la base.

Los datos obtenidos como resultado de este método incluyen el valor base, el tiempo necesario para alcanzar el valor base, el tiempo necesario para alcanzar la suma de los valores base y de comienzo, y la diferencia de estos dos períodos de tiempo (tiempo Base hasta Iniciación).

Con valores de medición cinética en aumento, el valor de iniciación se debe definir como un número positivo, mientras que con valores decrecientes el valor de iniciación se debe definir como un número negativo.

Mín./Máx./Área

Utilice esta pestaña para definir la evaluación de los valores mínimo y máximo contenidos en la curva y para definir el cálculo del área bajo la curva cinética:

Lista desplegable Datos de entrada	Seleccione los datos de entrada que desea procesar.
Casilla de verificación Valor mínimo de la media	Selecciónelo para abrir el campo de texto Puntos donde se escribe el número fijo de puntos para el valor mínimo. Entonces se usarán esos puntos para generar una curva uniforme y se determinará el valor más bajo contenido en la curva.



Casilla de verificación Valor máximo de la media	Selecciónelo para abrir el campo de texto Puntos donde se escribe el número fijo de puntos para el valor máximo. Entonces se usarán esos puntos para generar una curva uniforme y se determinará el valor más alto contenido en la curva.
Casilla de verificación Área	En los campos de texto suministrados se pueden indicar los tiempos de inicio y fin, a no ser que se desee analizar toda la cinética.
	 Botones de opción Tiempo/Puntos: si se selecciona Tiempo, se deben escribir los tiempos de inicio y fin en horas, minutos y segundos. si se selecciona Puntos, se debe escribir el número de ciclos cinéticos para indicar los puntos de inicio y de final del análisis.
	 Campo Inicio: aquí se debe escribir la hora de comienzo (en horas, minutos y segundos) o el número de ciclos.
	 Campo Final: aquí se debe escribir la hora de finalización (en horas, minutos y segundos) o el número de ciclos.

Las medias se toman a partir del número de puntos desde el primer valor de medición cinética y en pasos de uno hasta que se incluya en el cálculo de la media el último punto de medición cinética. El valor mínimo/máximo se determina a partir de estas medias.

Los datos resultantes de este método incluyen los valores mínimo/máximo y el período de tiempo desde la primera medición hasta los valores mínimo/máximo en segundos.

Ejemplo:

Con una cinética de 5 ciclos y un número fijo de 3 puntos, se toman los siguientes promedios:

promedio del 1.º, 2.º y 3.º valor cinético

promedio del 2.º, 3.º y 4.º valor cinético

promedio del 3.º, 4.º y 5.º valor cinético

El valor mínimo/máximo se determina a partir de estas 3 medias.

El área bajo la curva se calcula utilizando la siguiente fórmula:

$$A = \sum_{i=1}^{n-1} y_i * (x_{i+1} - x_i) + \frac{(y_{i+1} - y_i) * (x_{i+1} - x_i)}{2}$$

Pestaña Datos disponibles

En la pestaña Datos de salida disponibles se muestra la lista de resultados:

Campo El campo de datos muestra los valores seleccionados en Datos El campo de datos muestra los valores seleccionados en las pestañas anteriores del cuadro de diálogo Parámetros de cálculos cinéticos. Este campo de datos únicamente sirve para mostrar los datos de salida, no dispone de funciones de edición.



Pestaña Cinética de enzima

Utilice esta pestaña para establecer la evaluación de cinéticas de enzimas según el modelo de Michaelis y Menten:

Lista desplegable Datos de entrada	Seleccione los datos de entrada que desea procesar.
Casilla de verificación Calcular Km y Vmáx	Seleccione la casilla de verificación correspondiente para indicar si desea calcular Km y Vmáx.
Cuadro de grupo Tipo de cálculo	 El tipo de cálculo se puede seleccionar como: Hanes (concentración frente a concentración/datos de entrada) Eadie-Hofstee (datos de entrada/concentración frente a datos de entrada)
	 Lineweaver-Burk (1/datos de entrada frente a 1/concentración)

Los datos resultantes de este método incluyen Km y Vmáx para el gráfico de cinética de enzimas de cada grupo experimental.

A diferencia de los resultados de los otros cálculos cinéticos, estos resultados se muestran en el cuadro de diálogo **Gráfico: Cinética de enzima**.

Esta pestaña solamente estará disponible si en la placa hay estándares y se han definido transformaciones o cálculos de pendiente.

4.3.7 Transformaciones cinéticas: Añadir nuevas transformaciones cinéticas

En el cuadro de diálogo **Transformaciones cinéticas** se puede definir una fórmula de transformación que se utilizará para transformar datos de entrada cinéticos individualmente para cada pocillo.

En el cuadro de diálogo **Transformaciones cinéticas** se pueden realizar cálculos adicionales sobre los datos cinéticos de entrada.

Los elementos de la ventana son similares a los de la entrada **Transformación**. Para obtener información adicional, consulte el capítulo 4.3.5 Datos transformados: Añadir nueva transformación.

4.3.8 Concentraciones: Curva estándar

Utilice esta opción para configurar curvas estándar para pruebas cuantitativas.

El cuadro de diálogo de la curva estándar contiene varios ajustes relativos al tipo de análisis, los ejes y la visualización de la curva estándar. Los campos y elementos editables están agrupados en cinco pestañas diferentes.



Pestaña Datos

Esta pestaña establece algunos ajustes básicos como el origen de los datos de entrada.

Control de entrada numérica Grupo exp.	Si se van a realizar varias pruebas sobre una placa, se debe seleccionar el Grupo experimental . Si la placa contiene solamente una prueba, entonces no se mostrará el control de entrada numérica. Para cada grupo experimental, los datos de entrada se pueden definir individualmente.
Lista desplegable Datos de entrada	Seleccione los Datos de entrada que se utilizarán para la curva estándar. Seleccione datos de medición o los resultados de cualquier transformación disponible.
Botón de opción Estándares de diseño	Calcula la curva estándar a partir de los estándares que hay en el diseño.
Botón de opción Estándares en archivo ext.	Si el grupo experimental seleccionado no contiene ningún estándar, se puede cargar una curva estándar desde un archivo .std. Para seleccionar el archivo, haga clic en el botón Seleccionar .
Botón de opción Estándares en grupo exp.	Si el grupo experimental seleccionado no contiene ningún estándar, se puede usar una curva estándar de otro grupo experimental.
Botón de opción Sin gráfico de curva estándar	Si el grupo experimental seleccionado no contiene ningún estándar, se puede decidir no calcular ninguna concentración para este grupo experimental (en este caso se establece como predeterminado).
Botón Concentraciones adicionales	Haga clic en el botón Concentraciones adicionales para abrir el cuadro de diálogo Calcular concentraciones adicionales . Seleccione conjuntos adicionales de datos de entrada que se usarán para calcular concentraciones basadas en la curva estándar actual.
El cuadro de diálogo Calcular concentraciones adicionales contiene los siguientes elementos:	
l ista despleaable	Seleccione los datos de entrada para el cálculo de

Lista desplegable Datos de entrada	Seleccione los datos de entrada para el cálculo de concentraciones adicionales.
Lista Datos seleccionados	Esta lista contiene los nombres de los datos de entrada para el cálculo de concentraciones adicionales.
Botón Añadir	Haga clic en el botón Añadir para agregar a la lista Datos seleccionados los datos de entrada seleccionados actualmente en la lista desplegable Datos de entrada .
Botón Eliminar	Haga clic en el botón Eliminar para quitar los datos seleccionados de la lista Datos seleccionados .



Pestaña Tipo de análisis

Utilice esta pestaña para seleccionar el tipo de análisis. Consulte en la sección 11.4 Tipos de análisis de la curva estándar una descripción detallada de los tipos de análisis.

Botones de opción Tipo de análisis	Seleccione qué algoritmo de interpolación se debe emplear para calcular la curva estándar:
	Se enumeran y pueden seleccionarse todos los tipos de análisis proporcionados:
	1. Punto a punto
	2. Regresión lineal
	3. Regresión no lineal
	4. Spline cúbica
	5. Akima
	 Polinómico (incluye el campo para seleccionar el orden polinómico e incluye opciones de ponderación)
	 Cuatro parámetros (incluye el botón Más para determinar mínimo y máximo)
	 Algoritmo de Marquardt con cuatro parámetros (incluye el botón Más para definir las opciones de ponderación)
	 Algoritmo de Marquardt con cinco parámetros (incluye el botón Más para definir opciones de ponderación)
	10.LogitLog (incluye el botón Más para determinar mínimo y máximo)
Lista desplegable Escala de datos	El tipo de análisis se aplica a los valores escalados. Pueden seleccionarse los siguientes modos de escalado:
	 Lin(x)Lin(y): los ejes x e y usan escala lineal
	 Lin(x)Log(y): el eje x usa escala lineal y el eje y usa escala logarítmica
	 Lin(x)Log(y): el eje x usa escala logarítmica y el eje y usa escala lineal
	• Lin(x)Lin(y): los ejes x e y usan escala logarítmica



Aviso

Tenga en cuenta que si se selecciona la escala logarítmica para el eje x no será posible tener un valor de concentración de 0, ya que no es matemáticamente posible calcular el logaritmo de 0. Pero se puede utilizar un valor de, por ejemplo, 0,0000001.



Botón Más	Haga clic en el botón Más para ver las siguientes opciones:
	 Campo de selección numérico Curva estándar Si hay más de un grupo experimental para LogitLog y Cuatro parámetros se pueden seleccionar individualmente los valores Mín. y Máx. para cada curva.
	 Ponderación Seleccione Uso de pesos para utilizar uno de los siguientes métodos de ponderación:
	 Automática usando varianza Uso automático de pesos relativos Manual
	 Consulte la sección 11.4.14 Ponderación para ajuste de cuatro / cinco parámetros – ajuste polinomial / Marquardt Cuadro de diálogo Ponderación – Seleccione el botón Más si ha seleccionado "Automática usando varianza" => Se muestra el cuadro de diálogo de gestión de errores – Defina los factores de ponderación, cuál se debe utilizar si se producen uno de los dos casos de error (todas las réplicas son iguales o solo queda una réplica), que produce una varianza de 0 y por tanto conduce a un error de cálculo. Campos de texto Mín/Máx Mín/Máx permite al usuario definir el límite mínimo o máximo de la curva estándar para LogItLog y Cuatro parámetros.
Casilla de verificación Incluir (0,0)	Si se selecciona la opción Incluir , se añadirá (0,0) como punto estándar. Esto solo está permitido si se ha seleccionado Lineal como división para ambos ejes.
Casilla de verificación Extrapolación	La extrapolación aplica los cálculos de concentración a aquellos puntos base que se encuentran fuera del rango permitido, así como a aquellos que se encuentran dentro del rango. Por ejemplo, un factor de extrapolación de 3 fuerza un cálculo de concentración de valores entre mín- 2*(máx-mín) y máx+2*(máx-mín), donde mín y máx corresponden al valor de concentración mínimo y máximo de la curva estándar.
Campo numérico Factor de extrapolación	El campo de datos de factor de extrapolación solamente estará activo si se selecciona la casilla de verificación Extrapolación . Define los nuevos límites para el cálculo de concentración.

Pestaña Intercepciones

Utilice esta pestaña para calcular las concentraciones de los valores de Y seleccionados basándose en la curva estándar.

Control de entrada numérica Grupo exp.	Si hay disponibles varios grupos experimentales, se puede seleccionar el grupo apropiado.
Lista desplegable Entrada	En esta lista se muestran todos los datos disponibles para los cálculos.
Campo de lista Fórmula y nombre de intercepción	Primero se debe crear un nombre para la intercepción. Después se debe definir la fórmula de intercepción. Se puede introducir un valor numérico o una fórmula.
Operadores Entrada de fórmula	Los cuadros de selección permiten introducir fórmulas fácilmente, pues ofrecen gran variedad de funciones, operadores y variables.
	 Variables Todas las variables aceptadas por el software están disponibles utilizando el cuadro de lista.
	 Botón Haga clic en este botón para abrir el cuadro de diálogo Definir constantes.
	 Operadores Aquí se muestran todos los operadores aceptados por el software.
	 Funciones Aquí se muestran todas las funciones aceptadas por el software.

Ejemplo

IC50 para curva estándar:

Fórmula: (ST1_1+ST1_8)/2

valor del estándar más pequeño más valor del estándar más alto, dividido por 2



Pestaña Ejes

Utilice esta pestaña para definir el aspecto de los ejes.

Cuadro de grupo Eje X	En cada grupo están disponibles los siguientes elementos:
Cuadro de grupo Eje Y	Campo de texto Etiqueta : Introduzca una etiqueta de eje
	 Botón Color: Seleccione un color para todos los elementos del eje
	 Casilla de verificación Escala log.: Seleccione esta casilla para mostrar los ejes del gráfico logarítmicamente (no afecta al cálculo).
	 Botón de opción Autoseleccionar rango: El software determinará automáticamente el valor mínimo y máximo del eje
	 Botón de opción Rango: Se activan los campos numéricos Mín. y Máx Se pueden especificar el valor mínimo y máximo del eje.
	Casilla de verificación Retícula: Si se selecciona, se mostrará una retícula en el eje. El botón Color y la

para personalizar la retícula.

lista desplegable de estilos Línea se pueden usar

Pestaña Gráfico

Utilice esta pestaña para definir el aspecto del gráfico.

Cuadro de grupo Título	 Están disponibles los siguientes elementos: Campo de texto Etiqueta: Introduzca una etiqueta para el gráfico:
	 Botón Color: Seleccione un color para la etiqueta del gráfico.
Cuadro de grupo	Están disponibles los siguientes elementos:
Curvas	 Control de entrada numérica Grupo experimental: Seleccione el grupo experimental que personalizará (solo está disponible si se ha definido más de un grupo experimental).
	 Botón Color: Seleccione un color para la curva del grupo experimental seleccionado.
	 Lista desplegable Símbolo: Seleccione un símbolo para los puntos base del grupo experimental seleccionado.
	 Campo de texto Etiqueta: Introduzca una etiqueta para la curva del grupo experimental seleccionado.
	 Casillas de verificación Esconder curva: Si se selecciona, no se mostrará la curva.
	 Control de entrada numérica Ancho de línea: Seleccione la anchura de línea para el grupo experimental seleccionado.
Cuadro de grupo Fuente	Seleccione un tamaño de fuente Pequeño , Medio o Grande .



Cuadro de grupo Visualizar	Se puede seleccionar la visualización de los siguientes elementos:
	Leyenda : El color, el estilo del punto base y la etiqueta de la curva se muestran junto a información adicional (parámetros de tipo de análisis, coeficiente de correlación, etc.).
	Puntos base
	Intercepciones : Si se especifican intercepciones, se pueden mostrar y etiquetar en el gráfico.
	Barras de errores : Si los estándares se definen como réplicas, se muestra una barra que indica el rango más/menos la desviación estándar para cada punto

4.3.9 Transformaciones de concentraciones: Añadir nueva transformación de concentración

base.

En el cuadro de diálogo **Transformaciones de concentración** se puede definir una fórmula que se utilizará para transformar individualmente los datos de concentración de entrada para cada pocillo.

En el cuadro de diálogo **Transformaciones de concentración** se pueden realizar cálculos adicionales sobre los datos de concentración de entrada.

Los elementos de la ventana son similares a los de la entrada **Transformación**. Para obtener información adicional, consulte el capítulo 4.3.5 Datos transformados: Añadir nueva transformación.

4.3.10 Evaluar datos: Definición del corte

Utilice esta opción para clasificar datos sin procesar o datos calculados en límites de umbral. Los límites de umbral se pueden definir como valores numéricos fijos o como fórmulas.

Utilice la ventana **Definición del corte** para definir los rangos del corte y para asignar colores y nombres a resultados del corte.

La ventana Definición del corte contiene los siguientes elementos:

Lista desplegable Datos de entrada	Seleccione los datos de entrada que se van a utilizar en la evaluación.
Lista de selección Grupo exp.	Si la placa contiene más de una prueba, se debe seleccionar el grupo experimental correspondiente al que se va a aplicar el corte. Si la placa contiene solamente una prueba, solamente habrá disponible un grupo y ese campo no estará visible.

En la lista **Corte**, se pueden definir hasta diez rangos de corte. Se pueden escribir los límites para cada rango.

En la parte superior de la lista se debe escribir el valor límite superior. El siguiente valor inferior se indica a continuación, etcétera. Una flecha Bajo-Alto muestra el nivel. Los valores límites en sí pertenecen a los rangos superiores (condición mayor o igual).

Botones	Se puede asignar color a los niveles de resultados
Colores	individuales. Cuando se muestren los resultados, se
	aplicarán esos colores para representar los resultados
	cualitativos.

Campos de texto Etiqueta	Se debe asignar nombre a los indicadores de resultados para cada rango de valores, como por ejemplo: positivo (pos), negativo (neg), intermedio (?),
Campos de texto Límites	Los límites se pueden escribir como un valor constante o como una fórmula. Las fórmulas se pueden escribir directamente en un campo o se pueden usar campos Entrada de fórmula . Se puede definir un máximo de nueve límites.
Cuadro de grupo Entrada de fórmula	Las fórmulas que se van a escribir en el campo Límites se pueden componer mediante las variables, operadores y funciones siguientes. Para obtener detalles adicionales, consulte el capítulo 11.3 Escribir una fórmula.
	 Lista desplegable Variables Muestra todas las variables disponibles.
	 Botón Haga clic en este botón para abrir el cuadro de diálogo Definir constantes.
	 Lista desplegable Operadores Muestra todos los operadores disponibles.
	 Lista desplegable Funciones Permite seleccionar las funciones disponibles.
Casilla de verificación Prueba competitiva	Las pruebas competitivas asignan un resultado positivo a los valores bajos y un resultado negativo a los valores altos. En la evaluación, los límites irán de arriba abajo y se asignará el símbolo de los resultados correspondientes cuando un valor sea menor que o igual que el límite. En vista de ello, se invertirá la flecha Bajo-Alto .
Botón Seleccionar resultados de corte	Haga clic en el botón Seleccionar resultados de corte para abrir el cuadro de diálogo Seleccionar resultados de corte.

Seleccionar resultados de corte

Utilice esta ventana para seleccionar si se mostrará o no un resultado cualitativo para un cierto tipo de identificador.

La ventana **Seleccionar resultados de corte** contiene los siguientes elementos:

Estructura de árbol	Se muestran en una estructura de árbol todos los nombres
Selección de	de identificadores definidos, agrupados por tipo de
visualización de	identificador. Cada identificador va asociado a una casilla
cortes	de verificación.
	Al seleccionar o deseleccionar la casilla de verificación de un nombre de identificador se indica si se mostrarán o no los resultados cualitativos de los pocillos correspondientes. Después de la evaluación, solamente se mostrarán los resultados de corte para los tipos de identificadores seleccionados.



4.3.11 Evaluar datos: Validación QC

Las validaciones se utilizan para verificar la validez de una prueba. Si no se cumplen los criterios definidos, se mostrará un mensaje de error después de la ejecución de la medición y no se mostrará ningún resultado de corte (excepto si el usuario tiene el derecho **continuar evaluación en caso de error**. Consulte el capítulo 9.6 Derechos del usuario).

Una vez que se seleccionan los datos de entrada correctos, se deben escribir las fórmulas deseadas en el campo de lista **Condiciones de validación**. Un editor de fórmulas ayuda al usuario a elaborar las fórmulas matemáticas correspondientes.

Ejemplo:

Si los criterios de validación se definen como sigue:

NC1>0 y NC1<0,1

la evaluación comprobará si el control negativo se puede encontrar dentro del rango dado. Si no es así, se mostrará un mensaje de error.

El cuadro de diálogo **Definir validaciones del QC** contiene los siguientes elementos:

Lista desplegable Datos de entrada	Seleccione los datos correspondientes a los cuales debe aplicarse la validación. Por ejemplo: Datos sin procesar, etc.
Campo de selección Grupo exp.	Si la placa contiene más de una prueba, se debe seleccionar el grupo experimental correspondiente al que se va a aplicar la validación. Si la placa contiene solamente una prueba, solamente habrá disponible un grupo.
Campo de selección Grupo de validación	Use Grupos de validación para definir criterios de validación para el mismo grupo experimental con datos de entrada diferentes.
Campo de lista Condiciones de Validación	Esta lista unidimensional se rellenará con las fórmulas y las ecuaciones lógicas que definen las Condiciones de validación . Estas ecuaciones lógicas generarán un resultado lógico. El programa usará esa ecuación para examinar los datos de entrada y, si se cumplen los criterios de validación, devolverá el resultado VERDADERO. Si no se satisfacen los criterios, se mostrará el resultado FALSO y un mensaje de error.
Cuadro de grupo Entrada de fórmula	 Partes de las fórmulas que se escribirán en el campo Condiciones de validación se pueden seleccionar en las siguientes listas desplegables: Lista desplegable Variables Muestra todas las variables disponibles. Botón Haga clic en este botón para abrir el cuadro de diálogo Definir constantes. Lista desplegable Operadores Muestra todos los operadores disponibles. Lista desplegable Funciones A través de la paleta dada se pueden seleccionar las funciones disponibles. Para obtener detalles adicionales, consulte el capítulo 11.3 Escribir una fórmula .



QC de placa a placa

Utilice esta opción para definir una validación QC de largo plazo (consulte el capítulo 7.4.2 Menú de Barra de herramientas: Archivo/ QC de placa a placa):

Lista desplegable Datos de entrada	Seleccione en la lista desplegable los datos relevantes a los cuales se debe aplicar la validación.
Casillas de selección Control	Seleccione en la lista desplegable el tipo de control y luego escriba la media y la desviación estándar previstas.
	Para definir más de cuatro controles de QC se muestra una barra de desplazamiento junto a los controles del cuadro de diálogo.
Opciones de Zonas de trabajo para evaluar	Seleccione el filtro apropiado para definir qué zonas de trabajo deberían ser evaluadas.

4.3.12 Procesamiento de datos: Exportación de datos

Mediante este cuadro de diálogo se pueden seleccionar datos para exportarlos a archivos ASCII o Excel:

Lista Datos disponibles	Se muestra una lista de todos los datos disponibles, en función de la definición del método.
\leftarrow y \rightarrow (botones)	Los datos para exportar se pueden seleccionar simplemente con el método de arrastrar y colocar, o bien, haciendo clic en los datos de la ventana Datos disponibles y luego en la flecha que señala hacia la ventana Datos seleccionados . Para anular la selección de los datos, se puede seguir el procedimiento inverso.
Arriba y Abajo (botones)	Se puede seleccionar un elemento y pulsar el botón Arriba para subirlo o Abajo para bajarlo y así cambiar el orden de los datos seleccionados.
Lista Datos seleccionados	Esta lista muestra todos los datos seleccionados, transferidos desde la lista de campos Datos disponibles.
Botón Opciones al exportar	Haga clic en este botón para abrir el cuadro de diálogo Opciones de exportación .

La exportación real de los datos solamente se produce mediante procesamiento automatizado de datos (consulte 4.3.14 Procesamiento de datos: Procesamiento automatizado de datos).



Opciones al exportar

El cuadro de diálogo **Opciones al exportar** contiene los siguientes elementos:

Cuadro de grupo Dirección	El usuario puede definir si los datos de placa se extraen horizontalmente línea a línea o verticalmente columna a columna y si se escriben en ese orden en el archivo.	
	 Botón de opción Horizontal Los datos se recopilan en filas. 	
	 Botón de opción Vertical Los datos se recopilan en columnas. 	
Cuadro de grupo Resultado	Seleccione cómo exportar los datos, en forma de matriz o como una lista continua:	
	Botón de opción Matriz (anidada) Los datos de todos los conjuntos de datos seleccionados se organizan en una sola matriz. Esta matriz contiene las primeras columnas de todas las series de datos seguida por las segundas columnas de todas las series de datos.	
	 Botón de opción Matriz (separada) Los datos de cada conjunto de datos seleccionado se organizan en una matriz independiente. 	
	 Botón de opción Matriz (estilo XFluor) Los datos de cada conjunto de datos seleccionado se organizan en una matriz independiente junto con descripciones de pocillos y columnas similares a XFluor. 	
	 Botón de opción Tabla (pocillos en filas) Todos los datos de un pocillo se organizan en dirección horizontal empezando por los datos de la posición de pocillo A1. 	
	 Botón de opción Tabla (datos de pocillos en columnas) Todos los datos de un pocillo se organizan en dirección vertical empezando por los datos de la posición de pocillo A1. 	
	Si se selecciona la casilla de verificación Añadir marcas de tiempo cinéticas, se agregan las marcas de tiempo de las mediciones. Si se selecciona la casilla de verificación Añadir temperaturas, se agregan las temperaturas de las mediciones	
	Si se selecciona la casilla de verificación Insertar nombres datos , se muestra una lista con los nombres de los datos exportados encima del archivo ASCII o la hoja de cálculo Excel respectivos.	
Cuadro de grupo Añadir datos	Proporciona opciones de datos seleccionables. La información seleccionada se adjuntará al final de los datos y después se exportará con los datos seleccionados. Seleccione las casillas correspondientes:	



	Fecha/hora de medición
	Nombre de archivo del método
	Ruta de acceso del método
	Nombre de archivo de la zona de trabajo
	Ruta de acceso de la zona de trabajo
	 Valor(es) de longitud de onda del filtro
	Indicaciones al usuario
	Nombre del usuario actual
	Parámetros de medición
	 Información de placas múltiples
	 Registro de modificaciones de la zona de trabajo (solo magellan Tracker)
	 Firmas de la zona de trabajo (solo magellan Tracker)
Botón Establecer como predeterminado	Los ajustes se pueden registrar como valores predeterminados para su uso futuro.
Botón Restablecer valores predeterminados	Los ajustes se pueden restablecer a los valores predeterminados definidos previamente.

Ejemplo

Datos sin procesar, ciclo cinético 1, marca de tiempo 0 s

11 12 13

21 22 23

Datos sin procesar, ciclo cinético 2, marca de tiempo 33 s

81 82 83

91 92 93



Matriz (anidada), horizontal: 11 81 12 82 13 83 21 91 22 92 23 93	Matriz (separada), horizontal, con marcas de tiempo: 0 s 11 12 13 21 22 23 33 s 81 82 83 91 92 93	Matriz (estilo XFluor): <> 1 2 3 A 11 12 13 B 21 22 23 <> 1 2 3 A 81 82 83 B 81 82 83
Tabla (datos de pocillos en filas), horizontal, con marcas de tiempo: 0 s 33 s 11 81 12 82 13 83 21 91 22 92 23 93	Tabla (datos de pocillos en columnas), horizontal: 11 12 13 21 22 23 81 82 83 91 92 93	Tabla (datos de pocillos en columnas), vertical: 11 21 12 22 13 23 81 91 82 92 83 93

Exportar a archivo ASCII

El cuadro de diálogo **Exportar a archivo ASCII** contiene los siguientes elementos:

Cuadro de grupo Carácter decimal	 El carácter decimal puede ser un punto o una coma: Botón de opción Punto (x.xx) Botón de opción Coma (x,xx)
Cuadro de grupo Delimitador	Esta opción establece un delimitador para separar valores individuales de datos en las líneas de un archivo ASCII:
	 Botón de opción Tabulador Opción predeterminada, se utilizan tabuladores como delimitadores.
	 Botón de opción Otro símbolo Si lo selecciona, puede escribir un carácter delimitador definido por el usuario.
Cuadro de grupo Directorio	 Esta opción establece el directorio donde se guardará el archivo ASCII: Botón de opción Directorio predeterminado de exportación Botón de opción Utilizar directorio: Seleccione el directorio donde desea guardar el archivo ASCII
Cuadro de grupo Codificación	 Seleccione el código de exportación: ANSI: Genera caracteres compatibles con ANSI. Unicode: Para caracteres que no se pueden exportar con código ANSI (p. ej., chinos o cirílicos).


Botón Establecer como predeterminado	Estos ajustes se pueden registrar como predeterminados para su uso futuro.
Botón Restablecer valores predeterminados	Los ajustes se pueden restablecer a los valores predeterminados definidos previamente.



SE DEBE SELECCIONAR UN DELIMITADOR DE EXPORTACIÓN APROPIADO. EL DELIMITADOR DE EXPORTACIÓN Y EL CARÁCTER DECIMAL NO DEBERÍAN SER EL MISMO. EL TABULADOR ES EL DELIMITADOR PREDETERMINADO, PERO SE PUEDE SELECCIONAR OTRO SÍMBOLO.

Exportación a Excel

El cuadro de diálogo Exportar a Excel contiene los siguientes elementos:

Cuadro de grupo Objetivo	 Los botones de opción definen dónde colocar los datos transferidos dentro de Excel. Esta opción también es válida para la exportación automática a Excel. Botón de opción Nuevo libro de trabajo Los datos transferidos se guardarán en la primera hoja de cálculo de un nuevo libro de trabajo. Siempre se creará un nuevo archivo, independientemente de si Excel está abierto o no.
	 Botón de opción Nueva hoja de cálculo Los datos transferidos se insertarán en una nueva hoja de cálculo de un libro Excel abierto y activado. Si Excel no está abierto, si se selecciona esta opción, se abrirá Excel, se creará un nuevo libro y los datos solicitados se insertarán en la primera hoja de cálculo.
	 Botón de opción Insertar en hoja de cálculo en la celda y campo de texto la celda coordenadas de la celda El primer valor de los datos transferidos se colocará en la celda indicada (A1 es la celda predeterminada) de una hoja de cálculo Excel abierta y activa. Si se ha seleccionado esta opción y Excel no está abierto, se iniciará Excel, se creará un nuevo libro, y se insertarán los datos solicitados en la primera hoja de cálculo, en la celda definida.
	 Botón de opción Anexar a la hoja de cálculo actual Esta opción añadirá los datos a la hoja de cálculo actual. Si se ha seleccionado esta opción y Excel no está abierto, se iniciará Excel, se creará un nuevo libro y se insertarán los datos solicitados en la primera hoja de cálculo.



	 Utilice Insertar en plantilla para exportar los datos a una plantilla Excel predefinida. Haga clic en el botón para escribir el directorio y el nombre de la plantilla Excel. Se debe indicar la correspondiente posición de celda.
Botón Establecer como predeterminado	Estos ajustes se pueden registrar como predeterminados para su uso futuro.
Botón Restablecer valores predeterminados	Los ajustes se pueden restablecer a los valores predeterminados definidos previamente.

4.3.13 Procesamiento de datos: Informe impreso

El cuadro de diálogo **Informe impreso** contiene funciones de formateo para impresoras de copias impresas que permiten personalizar el contenido y el aspecto de la impresión.

Los ajustes de los informes están agrupados en cuatro pestañas, que representan distintos aspectos del formateo:

- Pestaña Selección de datos
- Pestaña Preparar página
- Pestaña Encabezado
- Pestaña Pie de página

Pestaña Selección de datos

La pestaña **Selección de datos** ofrece una lista de datos disponibles para imprimir.

Cuadro de grupo Imprimir como	Seleccione si los datos se imprimirán en forma de matriz o en forma de lista:		
	 Lista El formato de lista individual (horizontal, vertical, ordenada y alternada) se debe configurar en el cuadro de diálogo de propiedades (véase a continuación). 		
	 Matriz Al seleccionar matriz los datos se imprimen en una lista bidimensional (matriz). 		
Listas de origen y destino de los datos	 Para definir los datos impresos se usan dos listas: Lista Datos disponibles: Lista de todos los datos disponibles que se pueden imprimir. Se pueden agregar saltos de página y líneas divisorias para obtener el diseño deseado. Se puede hacer doble clic para seleccionar los elementos que se imprimirán. Se enumeran en la lista de datos seleccionados. 		

	 Lista Datos seleccionados: Lista de todos los datos seleccionados que se imprimirán. Las notas informativas también ayudan a dar una idea de cómo se imprimirán los datos. Por ejemplo, pueden indicar si los datos se mostrarán en forma de matriz, lista, recopilación de datos o como un gráfico. Las entradas que se muestran bajo un encabezado se incorporarán al formato de lista o matriz correspondiente, lo cual ofrece la posibilidad de comparar distintos conjuntos de valores, como por ejemplo valores de medición y resultados.
Botón Anexar	Haga clic en este botón para transferir los datos resaltados en la ventana de datos Disponibles a la ventana de datos Seleccionados . También puede utilizar arrastrar y colocar .
Botón Insertar	Use esta opción para mostrar dos secciones de datos con formato similar dentro de la misma área. En primer lugar, deben seleccionarse los datos de la lista Datos disponibles y después los datos de la lista Datos seleccionados . Al hacer clic en el botón Insertar , los primeros se incluirán dentro de los últimos y entonces ocuparán la misma área en la copia impresa que su elemento de origen. Por ejemplo, si los valores de desviación estándar se incluyen en el elemento de valor de medición, ambos se mostrarán dentro de la celda correspondiente de una matriz. En este caso, los valores de medición servirán como elemento de origen y serán, por lo tanto, los que se muestran primero.
Botón Quitar	Se pueden borrar los elementos seleccionados dentro de la lista Datos seleccionados .
Botones Arriba y Abajo	Use estos botones para reorganizar la secuencia de listas, gráficos, matrices, en la copia impresa o para reorganizar elementos dentro de una matriz o una lista.
Botón Propiedades	Se activa cuando se selecciona una lista. Se puede seleccionar el estilo y el contenido de la lista.
Cuadro de grupo Estilo	Botón de opción Lista vertical Imprime los datos en forma de lista vertical (orden A1, B1, etc.).
	Botón de opción Lista vertical alternada El orden de los datos se alterna para cada fila o columna (orden A1, B1, H1, H2, G2, A2)
	Botón de opción Lista horizontal Imprime los datos en forma de lista horizontal (orden A1, A2, etc.).
	Botón de opción Lista horizontal alternada El orden de los datos se alterna para cada fila o columna (orden A1, A2,A12, B12, B11,B1).



Botón de opción Lista ordenada Esta opción ordenará las entradas de todas las columnas conforme a las entradas de la primera columna. Las entradas alfanuméricas de la primera columna se dividen en dos partes: la primera solo contiene letras, la segunda solo números. Las entradas primero se ordenan por orden alfabético. Las entradas con el mismo nombre después se ordenan numéricamente. Ejemplo: Puede utilizar esta opción con los ID de muestras o con el diseño, restringido a un grupo experimental. La ordenación de los valores numéricos solo funcionará si todas las entradas son números
enteros.
Botón Lista(s) para muestras Permite la impresión de listas individuales por muestra. La separación para las tablas simples se realiza por los ID de las muestras (si están disponibles) o por nombres

ID de las muestras (si están disponibles) o por nombres de identificadores del diseño de placa. Se necesitan los ID de muestras si el diseño contiene distintos grupos experimentales que representan distintos parámetros para las mismas muestras. Se puede combinar el estilo de **Lista(s) para muestras** con otros **parámetros** como Identificadores, Grupos exp. y resultados de corte.

tilo omolou				Limitaciones	ОК
Layo Raw Blar Cuta 1 A1 A2 Layo Raw Blar	out data the reduction off results 2 BL1_1 0.1 BL1_2 0.1 out data the coults	3 4 101~ -0.000 102~ 0.000	 Lista vertical Lista vertical alternada Lista horizontal Lista horizontal alternada Lista ordenada Lista(s) para muestras 	Identificador Structure Structure Structure Structure Structure Structure Structure Structure Structure Structure Structure Structure Structure Structure Structure Structure Structure Structure Structure Structure Structure Structure Structure Structure Structure Structure Structure Structure Structure Structure Structure Structure Structure Structure Structure Structure Structure Structure Structure Structure Structure Structure Structure Structure Structure Structure Structure Structure Structure Structure Structure Structure Structure Structure Structure Structure Structure Structure Structure Structure Structure Structure Structure Structure Structure Structure Structure Structure Structure Structure Structure Structure Structure Structure Structure Structure Structure Structure Structure Structure Structure Structure Structure Structure Structure Structure Structure Structure Structure Structure Structure Structure Structure Structure Structure Structure Structure Structure Structure Structure Structure Structure Structure Structure Structure Structure Structure Structure Structure Structure Structure Structure Structure Structure Structure Structure Structure Structure Structure Structure Structure Structure Structure Structure Structure Structure Structure Structure Structure Structure Structure Structure Structure Structure Structure Structure Structure Structure Structure Structure Structure Structure Structure Structure Structure Structure Structure Structure Structure Structure Structure Structure Structure Structure Structure Structure Structure Structure Structure Structure Structure Structure Structure Structure Structure Structure Structure Structure Structure Structure Structure Structure Structure Structure Structure Structure Structure Structure Structure Structure Structure Structure Structure Structure Structu	Cancela Ayuda
1	2	3			
A3 A4 B3 B4	SM1_2_1 SM1_2_3 SM1_2_2 SM1_2_4	0 103~ 0.104~ 0 203~ 0 204~		Resultados de corte Todos los resultados de corte Solamente	

Cuadro de grupo Limitaciones Seleccione una de estas opciones:

Identificadores: seleccione o desmarque qué identificadores se deben mostrar en la lista.

Grupo(s) exp.: seleccione entre todos los grupos experimentales o introduzca el número de grupo experimental que desea mostrar en la lista.

Resultados de corte: seleccione entre todos los resultados de corte o seleccione en la lista qué resultados deben mostrarse en la lista.



Pestaña Preparar página

Cuadros de grupo Mostrar encabezado y pie de página	 Casilla de verificación Imprimir encabezado Si se selecciona, se imprimirá el encabezado Casilla de verificación Imprimir pie de página Si se selecciona, se imprimirá el pie de página Botones de opción de Encabezado Todas las páginas (imprime el encabezado en cada página) Solo la primera página (imprime el encabezado solo en la primera página) Excepto la primera página (imprime el encabezado en todas las páginas excepto la primera) Botones de opción de Pie de página Todas las páginas (imprime el pie de página en cada página) Excepto la primera página (imprime el pie de página en cada página) Botones de opción de Pie de página solo en la primera página (imprime el pie de página en cada página) Solo la primera página (imprime el pie de página solo en la primera página (imprime el pie de página solo en la primera página (imprime el pie de página solo en la primera página)
Grupo imprimir/página	 Botón de opción Color Si lo admite la impresora, la impresión será en color. Los datos específicos del pocillo se imprimen en el color del diseño (por ejemplo: los datos sin procesar de la muestra se imprimen en azul). Los resultados de corte se imprimen en el color del rango (por ejemplo: los pocillos "pos" se imprimen en rojo). El gráfico imprime tal como se muestra. Botón de opción Blanco y negro La impresión será en blanco y negro.
Cuadro de grupo Márgenes	 En la lista desplegable Unidad se pueden seleccionar centímetros o pulgadas. Hay 4 campos numéricos que definen los márgenes: Campo de texto Izquierda: introduzca el valor del margen izquierdo. Campo de texto Derecha: introduzca el valor del margen derecho. Campo de texto Sobre el texto: introduzca el valor del margen superior. Campo de texto Bajo el texto: introduzca el valor del margen inferior.

En la pestaña **Preparar página** se define la configuración general:



Pestaña Encabezado

En la pestaña **Encabezado** se define el contenido del encabezado del documento:

Selección de datos	Configuración Encabe	zado Pie de pág	gina			
0	oción:	Ajuste:			Número de fila:	
] Paginación] Fecha] Tiempo] Nombre del archivo] Nombre del programa] Comentario	 izquierda izquierda izquierda izquierda izquierda izquierda izquierda 	Centro centro centro centro centro	derecha derecha derecha derecha derecha derecha derecha derecha	3 \$\overline\$ 2 \$\overline\$ 3 \$\overline\$ 1 \$\overline\$ 1 \$\overline\$ 1 \$\overline\$	
	Separadores	✓ Inferior				
Casillas de verificación	Use cas	e Opción illas de ve	para se erificaci	elecciona	ar cualquiera de las s que se impriman en	iguientes
Opción	enc • 1	abezado: Paginació	: òn			
	• 1	Fecha (ad	ctual)			
	• -	⊣ora Nombre c	lel archi	vo		
	•	Nombre c	del prog	rama		
	. (Comenta	rio			
	El c sele	ampo de ecciona la	texto C a casilla	omenta de verif	rrio estará activo si s icación Comentario.	е
Botones de o	pción Sele	eccione u	ına de e	stas opo	ciones:	
Ajuste	• 1	zquierda				
	• (Centro				
	•	Derecha				
₋ista Número de f i	Núr Ia esta	nero de f ablecer la	fila es u fila par	in campo a cada c	o numérico que pern opción.	nite



Casillas de	Las casillas de verificación de separadores definen dónde
verificación	aparecerán las líneas delimitadoras. Se pueden colocar
Separadores	encima, debajo o encima y debajo del encabezado.
-	Cosillo de verificación Superior

- Casilla de verificación Superior: La línea de separación se muestra encima del encabezado.
- Casilla de verificación Inferior: La línea de separación se muestra debajo del encabezado.

Pestaña Pie de página

En la pestaña **Pie de página** se define el contenido del pie de página del documento. La pestaña **Pie de página** contiene los mismos elementos que la pestaña **Encabezado**. Para más detalles sobre la pestaña Pie de página, consulte la pestaña Encabezado.

La pestaña de pie de página desaparecerá si no se selecciona **Imprimir pie de página** en la pestaña **Preparar página**.

4.3.14 Procesamiento de datos: Procesamiento automatizado de datos

Después de una medición se pueden iniciar automáticamente algunas acciones.

Seleccione esas acciones y se ejecutarán cronológicamente en el siguiente orden:

- Si está seleccionada la casilla de verificación Cargar lista de IDs de muestras, se cargará automáticamente una lista de IDs de muestras.
- Si está seleccionada la casilla de verificación Guardar zona de trabajo (siempre está seleccionada en magellan Tracker), se guardará automáticamente la zona de trabajo creada.
- Si está seleccionada la casilla de verificación Exportar a archivo ASCII, los datos seleccionados para exportar se exportarán automáticamente a un archivo ASCII.
- Si está activada la casilla de verificación Exportar a archivo ASTM, después de ejecutar el método los datos seleccionados para exportar se exportarán automáticamente a formato ASTM para poderlos procesar en sistemas compatibles con LIS (Laboratory Information Systems).
- Si está activada la casilla de verificación Exportar a rastreo de muestras, después de ejecutar el método, los datos seleccionados para exportar se exportarán automáticamente al sistema de rastreo de muestras. Para obtener más información, consulte el manual correspondiente de rastreo de muestras.
- Si está seleccionada la casilla de verificación Exportar a Excel, los datos seleccionados para exportar se exportarán automáticamente a Excel.
- Si está seleccionada la casilla de verificación Imprimir, los datos seleccionados para el informe impreso se imprimirán automáticamente en la impresora predeterminada.
- Si está seleccionada la casilla de verificación Ver resultados después de la medición, se mostrarán para su examen los datos de mediciones y los datos calculados.



Más – Cargar lista de IDs de muestras

El cuadro de diálogo **Más (Cargar lista de IDs de muestras)** contiene los siguientes elementos:

Cuadro de grupo Nombre de archivo	Seleccione el nombre de archivo de la lista de IDs de muestras que desea cargar:
	 Botón de opción Mismo nombre de archivo que la zona de trabajo
	 Botón de opción Nombre de archivo: escriba un nombre de archivo en el campo de texto.
Cuadro de grupo Directorio	Seleccione el directorio donde está almacenada la lista de IDs de muestras:
	 Botón de opción Directorio predeterminado de listas de IDs de muestras
	 Botón de opción Utilizar directorio: Seleccione el directorio que desee.
Cuadro de grupo Formato	Seleccione el formato de lista de IDs de muestras que desea cargar (consulte los detalles en el capítulo 5.2.3 Importar lista de IDs de muestras):
	 Botón de opción Detección automática desde la lista de formatos admitidos
	 Botón de opción Formato de archivo personalizado: Haga clic en el botón Formato de archivo para abrir el cuadro de diálogo Formato especificado por el usuario.



Si se carga automáticamente una lista de IDs de muestras llena solamente de modo parcial, se eliminarán automáticamente del diseño todas las muestras no usadas. Si en el diseño se ha definido un control final (p. ej., en el pocillo H12), ese control se moverá automáticamente al pocillo a continuación del último pocillo de muestra utilizado.



Aviso

Las listas de IDs de muestras cargadas automáticamente solamente deben contener IDs para muestras y no deben contener IDs para no muestras (control, estándares, etc.).



Más – Nombre de Zona de trabajo

Al seleccionar **Guardar zona de trabajo – Más**, se muestra el cuadro de diálogo **Nombre de zona de trabajo** con los siguientes elementos:

Cuadro de grupo Formato	 Se debe seleccionar uno de los botones de opción disponibles para definir el nombre de la zona de trabajo predeterminada. Ese nombre puede constar de la fecha, un contador o un cierto conjunto de letras. Otra posibilidad es combinar la fecha con una serie de letras. Los contadores que se utilizan junto con la fecha se restablecen a diario; de lo contrario, el contador se incrementa con cada medición: Botón de opción Fecha (DDMMAAAA) + contador Botón de opción Fecha (AAAAMMDD) + contador Botón de opción Contador total Botón de opción Usar código de barras disponible:
Campo de texto Prefijo	El texto escrito se añade al comienzo del nombre de la zona de trabajo definido más arriba.
Campo de texto Ejemplo	El campo de texto Ejemplo muestra un ejemplo de la selección actual.
Botón E stablecer como predeterminado	Estos ajustes se pueden registrar como predeterminados para su uso futuro.
Botón Restablecer valores predeterminados	Los ajustes se pueden restablecer a los valores predeterminados definidos previamente.

Más – Exportar a LIS

El cuadro de diálogo Más (Exportar a LIS) contiene los siguientes elementos:

Cuadro de grupo	Defina los delimitadores ASTM:		
Definición de delimitador ASTM	 Delimitador de campo: escriba un signo separador de campos (por omisión) 		
	 Repetir delimitador: escriba el signo para repetir (\ por omisión) 		
	• Delimitador de componente : escriba un signo para separar componentes (^ por omisión).		
	 Carácter de escape: escriba el signo de escape (& por omisión). 		
Cuadro de grupo ID del remitente	Defina la configuración de ID del remitente que debe mostrar el encabezado de los archivos ASTM:		
	 Campo de texto ID del remitente: de modo predeterminado está vacío. Si se define, p. ej., SUNRISE, el nombre se muestra en el encabezado del archivo ASTM (p. ej. H \^& SUNRISE^Magellan V7.2^Tracker). 		
	 Anexar nombre de apl.: Si lo selecciona, en el encabezado del archivo ASTM se exportará el nombre de la aplicación (p. ej., Magellan V7.2[^]Tracker). 		



	 Anexar versión de apl.: Si lo selecciona, en el encabezado del archivo ASTM se exportará el número de versión de Magellan (p. ej., Standard o Tracker).
Cuadro de grupo Nombres de ensayos LIS	Seleccione el nombre del ensayo LIS que se mostrará en el registro de pedido de prueba y el registro de resultado del archivo ASTM:
	 Usar nombre del método: nombre del método que se utilizó para la medición.
	 Campo de texto Usar nombre: defina un nuevo nombre para mostrar.
	Ejemplo: ASTM_Test
	 Registro de pedido de prueba: O 1 ID_muestra ^^^nombre de método Fecha y hora de medición
	 Registro resultante: R 1 ^^^nombre_método^resultados
Cuadro de grupo	Seleccione el código de exportación:
Codificación	ANSI: Genera caracteres compatibles con ANSI.
	Unicode: Para caracteres que no se pueden exportar con código ANSI (p. ej., chinos o cirílicos).
Cuadro de grupo Directorio	 Defina el directorio para la exportación ASTM: Exportar directorio predeterminado: la exportación se realiza a la carpeta definida en Otras opciones.
	 Campo de texto Usar directorio: defina un nuevo directorio de exportación.
Botón Establecer como predeterminado	 Al usar esta opción, se pueden registrar los ajustes como predeterminados para su uso futuro.
Botón Restablecer valores predeterminados	 Al usar esta opción, se pueden restablecer los ajustes a los valores predeterminados definidos previamente.

Exportación ASTM – Archivo de ejemplo

Cada archivo de exportación ASTM (.txt) consta de los siguientes componentes:

- Registro de encabezado de mensaje: H|\^&|||SUNRISE^Magellan V7.2^Tracker
- Registro de datos de pacientes (P): P|1||ID_muestra 1 para el paciente 1
- Registro de solicitud de prueba (O):
 O|1|ID_muestra||^^^nombre de método||Fecha y hora de medición
- Registro resultante (R): R|1|^^^ nombre_método^resultados
- Registro de fin de mensaje: L|1|N



Ejemplo:

Configuración de exportación ASTM:

Definición de delimitador ASTM	ID del remitente		OK
Delimitador de campo: Image: Carácter de escape: Image: Carácter de escape: Carácter de escape: Image: Carácter de escape:	SUNRISE Anexar nomb aplicación Anexar versi aplicación	ore de ón de	Cancelar
Nombre LIS de ensayo		Codificación	
Directorio O Directorio predeterminado de	exportación		Establecer como predeterminado
Usar directorio: C:\Users\Public\Documents\	Tecan Magellan Jasc		Restablecer valores

Archivo ASTM correspondiente:

H|\^&|||SUNRISE^Magellan V7.2^Tracker P|1|| muestra1 O|1|muestra1||^^Prueba_ASTM||20180315091504 R|1|^^Prueba_ASTMt^Datos sin procesar^A1|0.101~|DO| P|2||muestra2 O|1|muestra2||^Prueba_ASTM||20180315091504 R|1|^Prueba_ASTM^Datos sin procesar^B1|0.201~|DO| P|3||muestra3 O|1|muestra3||^Prueba_ASTM||20180315091504 R|1|^Prueba_ASTM^Datos sin procesar^C1|0.301~|DO|



Más – Mostrar resultados

El cuadro de diálogo **Más (Mostrar resultados)** contiene los siguientes elementos:

Cuadro de grupo Vista de placa	Seleccione los datos que desea mostrar en la vista de placa:		
	 Lista desplegable Primera línea en pocillo (fija): Seleccione un nombre entre los nombres de datos disponibles. 		
	 Lista desplegable Segunda línea en pocillo (fija): Seleccione un nombre entre los nombres de datos disponibles. 		
	 Lista desplegable Tercera línea en pocillo (fija): Seleccione un nombre entre los nombres de datos disponibles. En contraste con las dos primeras líneas, esta selección solamente se aplica la primera vez. Siempre que se seleccionen en la barra de control otros datos para ser vistos en la vista de placa, esos datos se mostrarán en la 3ª línea. Las dos primeras líneas permanecen sin alterar. 		
	usar , que hace que la vista de placa resulte más legible.		
Cuadro de grupo Cuadro de texto	Seleccione un nombre entre los nombres de datos disponibles. Esta selección solamente se aplica la primera vez. Siempre que se seleccionen en la barra de control otros datos para ser vistos en el cuadro de texto, esos datos se mostrarán en el cuadro de texto.		

4.3.15 Otras opciones: Indicaciones al usuario

En la barra de control, expanda **Otras opciones** y seleccione **Indicaciones al usuario**.

En cada campo **Indicación** de la lista se puede introducir una palabra clave. Después se mostrará esa palabra clave antes de la medición. Se puede introducir un texto, como por ejemplo el número de lote o el nombre completo del analista.

Al seleccionar **Req.** (obligatorio), una medición solo se puede iniciar después de recibir una entrada (se debe añadir una información textual al campo **Respuesta a indicación**).

Las indicaciones también se mostrarán en la lista de datos disponibles para imprimir.

4.3.16 Otras opciones: Formato de números

En la barra de control, expanda **Otras opciones** y seleccione **Formato de números**.

El cuadro de diálogo **Formato de números** se usa para configurar el formato de los números al visualizar valores. Además, se puede seleccionar el formato científico para valores fuera de rango.

El ajuste predeterminado es 5 dígitos significativos. Al seleccionar **Definido por** el usuario se puede configurar un formato de número de 1 - 10 posiciones para enteros y 0 - 7 posiciones para decimales.

Si está seleccionado el formato de números definido por el usuario, es preciso configurar los dos campos de datos numéricos.



Además, se puede usar la notación científica (con exponente) o no científica. Se muestra un ejemplo del formato de la selección actual.

Haga clic en **Establecer como predeterminado** o en **Restablecer valores predeterminados** como corresponda para guardar la definición introducida.

4.3.17 Otras opciones: Notas de método

En la barra de control, expanda **Otras opciones** y seleccione **Notas de método**.

El campo **Notas de método** permite al usuario agregar al nombre de archivo algunos comentarios que se mostrarán en la copia impresa del método.

Se puede aplicar formato sencillo como texto en negrita, cursiva, subrayado y colores. Durante la impresión se puede omitir el formato.

4.4 Guardar métodos

Haga clic en **Finalizar** para guardar el método mediante la ventana **Guardar como**, que contiene los siguientes elementos:

Campo Nombre de archivo	Escriba o modifique el nombre de archivo del método.
Campo Comentarios del archivo	Añada algunos comentarios al nombre de archivo, que se mostrarán por ejemplo en los cuadros de diálogo Abrir archivo .
Campo Comentarios del registro de modificaciones	Añada algún comentario, que se guardará en el registro de modificaciones (solamente disponible para magellan Tracker).
Botón Organizar favoritos	Haga clic en este botón para abrir el cuadro de diálogo Organizar favoritos . Para conocer más detalles, consulte el capítulo 6.4 Activar favorito.
Casilla de verificación Ejecutar este método ahora	Seleccione esta casilla de verificación para que se inicie el método inmediatamente después de terminar este asistente.

Ese documento modificado se puede guardar con su nombre de archivo actual o con un nombre diferente.

Haga clic en **Guardar y finalizar** para guardar el método y cerrar el asistente **Crear o editar método**.

Botones adicionales de magellan Tracker:

Botón Registro de modificaciones	Haga clic en este botón para abrir el cuadro de diálogo Registro de modificaciones donde se muestran el usuario, la fecha, la acción ejecutada y un comentario.		
	Se puede hacer clic en los botones apropiados para ver una presentación preliminar de la zona de trabajo, la lista de métodos e IDs de muestras de las versiones anteriores y actual.		
	Además se puede guardar con otro nombre una versión anterior del método.		



Registro de modificaciones					
Usuario 1 Administrator (Admin) 2 Administrator (Admin)	Fecha 2009-07-21 10:29:40 2009-07-21 11:49:51	Acción Created Modified	Comenta	rio	
Presentación preliminar			Aceptar	Guardar versión prev	/ia como Ayuda
Botón Firma	Haga clic en e Firma , donde comentario. Este botón sol firmado el arch Para obtener r	ste botón p se muestra amente est nivo mediar nás detalle	ara abrir n: usuari á dispon nte el asis s, consul	el cuadro de o o, fecha, accio ible cuando se stente Anexar te el capítulo	diálogo ón y è haya ∲ firma .

8 Asistente Anexar firma.CampoCuando que escribe en una contraseña para un método,
ese método quedará protegido por contraseña y no se
podrá modificar hasta que se escriba la contraseña
correcta. Consulte la descripción detallada a
continuación.

4.4.1 Protección de métodos mediante contraseña

Proteger un método con contraseña

Los usuarios que tengan derecho para crear y para modificar métodos pueden asignarles una contraseña para protegerlos.

De manera predeterminada, los **Administradores** y los **Especialistas en aplicaciones** tienen los derechos necesarios.

Al crear un nuevo método, se puede definir la protección por contraseña en la pestaña **Guardar como** del asistente **Crear o editar método**. Añada el nombre del método, escriba la contraseña en el campo de edición **Contraseña de método** (esquina inferior derecha del cuadro de diálogo) y haga clic en **Finalizar**.

Confirme la contraseña en el cuadro de diálogo **Contraseña de método** que se muestra y haga clic en **Aceptar**. El método se guardará con la protección por contraseña.

Los métodos ya guardados pero que no están protegidos por contraseña se pueden abrir para protegerlos. Para ello, ábralos y defina una contraseña en la pestaña **Guardar como**, tal y como se ha descrito anteriormente.

Editar métodos protegidos con contraseña

Para abrir un método protegido por contraseña, escriba la contraseña de ese método. El método se abrirá en **modo Edición**. Podrá modificar y guardar el método.



Si se selecciona la opción **Solo lectura**, el método se abrirá en modo de solo lectura. En este caso también es posible modificar el método, pero no se podrán guardar los cambios.

Cambiar la contraseña de un método protegido con contraseña

Para cambiar la contraseña de un método protegido con contraseña, abra el método en **modo Edición** (es necesario escribir la contraseña) y escriba una nueva contraseña en el campo de edición **Contraseña de método** de la pestaña **Guardar como**. Haga clic en **Finalizar**.

Confirme la contraseña en el siguiente cuadro de diálogo **Contraseña de método** y haga clic en **Aceptar**.

4.5 Métodos multiplaca

Cuando se encuentra conectado un instrumento INFINITE F50 se pueden definir métodos multiplaca.

Durante la definición del método, en el editor de parámetros de medición se puede configurar el número de placas que se medirán.



Cuando sea preciso medir dos o más placas, también hay disponible en la ventana **Definir evaluación** un contador para cambiar entre placas.





Todos los parámetros adicionales se pueden definir como ya se ha visto para los métodos de placas simples. Al iniciar un método multiplaca se muestra el siguiente cuadro de diálogo:

jecutar medici	ión			
Medición de	placa	de 3		
Lista de IDs	de muestra:			Insertar
-				
	Parar	PAUSA	FIECUTAR	

Lista de IDs de muestra	Puede seleccionarse una lista de IDs de muestras para cada placa.
Botón PARAR	Se detiene la medición. Se muestran en la zona de trabajo creada los datos de las placas ya medidas y ya no es posible agregar mediciones de placas adicionales.
Botón PAUSA	La medición se detiene temporalmente. En la zona de trabajo creada se muestran los datos de las placas ya medidas, pero es posible agregar mediciones de placas adicionales.
Botón EJECUTAR	Se inicia la medición de la siguiente placa.
Introducir/Extraer placa	La placa se introduce o se extrae del instrumento.

Al ejecutar un método multiplaca se permite detener temporalmente la medición entre dos placas, guardar los resultados obtenidos y proseguir la medición para las placas restantes más adelante. Si se selecciona **Autoguardar zona de trabajo**, la zona de trabajo se guarda después de cada medición simple.

5. Asistente Crear o editar lista de IDs de muestras

5.1 Introducción

El asistente para **Crear o editar lista de IDs de muestras** se usa para crear nuevas listas de muestras o para modificar las listas de muestras existentes.

Resumen del flujo de trabajo

Arranque al asistente **Crear o editar lista de IDs de muestras**. Después de una página bienvenida, se muestra la página **Selección de archivo**. Seleccione una de estas opciones: **Crear nueva** lista de IDs de muestras o **Editar** una lista de IDs de muestras existente y guardar la lista de IDs a través de la página **Guardar como**.

5.2 Crear o editar lista de IDs de muestras

La ventana Selección de archivos contiene los siguientes elementos:

Botón de opción a) Crear nueva	Permite crear una nueva lista de IDs de muestras.	
Botón de opción b) Editar	Permite modificar las listas de IDs de muestras existentes.	
Cuadro combinado Mostrar	 Los archivos se muestran conforme a la selección: Todos los archivos Mis archivos: esta opción está disponible si está activada la administración de usuarios (siempre está activada en magellan Tracker). 	
Lista Nombre de archivo	Se debe seleccionar la lista de IDs de muestras que se va a editar. El campo de Notas junto a cada nombre de archivo contiene (si se ha rellenado) una breve descripción de la lista de IDs de muestras.	
	Se mostrarán todas las listas de IDs de muestras disponibles en el directorio de listas de IDs de muestras estándar que satisfagan los criterios de filtro seleccionados en el cuadro combinado Mostrar.	
Botón Presentación preliminar	Muestra una vista previa de la lista de IDs de muestras seleccionada y permite iniciar su impresión.	

5.2.1 Crear nueva lista de IDs de muestras

En la ventana **Seleccionar tipo de placa** se puede seleccionar el número de IDs de muestras por pocillo y el tipo de placa utilizado. La opción **Importar de lista externa de IDs de muestras** permite examinar distintos tipos de archivos de IDs de muestras e importarlos a **magellan**.



Aviso Algunas listas de IDs de muestras pueden contener información adicional tal como concentraciones y diluciones lo que podría afectar a la evaluación de los datos de medición.



Seleccione el tipo de placa Tipo de placa Tipo de placa desde método Qualitative_ELISA_Sunrise.mth Tipos de placa estándares 96 placa de pocillos Tipo de placa desde archivo pdl CLI961b.pdf O limportar desde lista externa de ID de muestras Importar	Configuración IDs de muestra por pocillo O Una Dos Irres Impresión Tabla horiz Tabla vert Matriz
Ayuda Cancelar <<< Atrás	ESCRIBIR IDs DE MUESTRAS

El cuadro de diálogo **Seleccionar tipo de placa** contiene los siguientes elementos:

Co	nfic	lura	ción	•
00	my	juia	CIUI	1

J	
Botones de opción IDs de muestras por pocillo	Escriba en el número de IDs de muestras requerido por pocillo (se puede indicar un máximo de tres).
Botones de opción Impresión	Defina el diseño para imprimir la lista de IDs de muestras.
Botones de opción Tipo de placa	 Cuatro botones de opción definen el tipo de placa que se puede seleccionar: Tipo de placa desde método Seleccione un método para cargar el diseño de placa. Haga clic en Cargar para examinar los archivos de métodos existentes. Se mostrará el cuadro de diálogo Abrir de Windows[™]. Seleccione el archivo que desee y haga clic en Abrir.
	 Tipos de placa estándares Para definir el formato de placa, se pueden seleccionar tipos placas de geometría estándar. Cuando se activa, en una lista desplegable asociada se puede seleccionar un tipo de placa de geometría estándar.
	• Tipo de placa desde archivo pdf Para definir el formato de placa se puede utilizar un archivo .pdf predefinido (instrumentos SUNRISE) o un archivo .pdfx (instrumentos INFINITE F50).
	Importar de lista externa de IDs de muestras Se puede seleccionar un tipo de archivo externo para importarlo como lista de IDs de muestras. Consulte el capítulo 5.2.3 Importar lista de IDs de muestras.



El botón **Escribir IDs de muestras** solamente estará habilitado si se han definido los ajustes requeridos (se ha cargado un método, se ha seleccionado un tipo de placa estándar o el archivo pdf).

5.2.2 Importar o modificar lista de IDs de muestras

Después de seleccionar el tipo de placa, se muestra la ventana de placa Importar o modificar lista de IDs de muestras:

- Barra de herramientas con botones de Configuración, Edición, Llenado automático, Eliminar, Deshacer/Rehacer, Formato de impresión, Presentación preliminar y Zoom
- · Vista de placa con un diseño esquemático de una microplaca
- Lista de IDs de muestras (cuando se haya seleccionado Tipo de placa desde método)
- Cuadro de diálogo para insertar IDs de muestras (cuando se haya seleccionado Tipo de placa estándar)

Si se va a crear una nueva lista de IDs de muestras, no se asignan IDs de muestras a los pocillos. Se puede hacer clic en un pocillo para seleccionarlo.

Es preciso escribir todos los IDs de muestras requeridos de modo manual o mediante los botones de la barra de herramientas, o bien se pueden leer con el escáner de códigos de barras portátil para IDs de muestras. Los IDs de muestras se pueden definir y editar en el tipo de microplaca definido. Se pueden introducir un máximo de tres IDs de muestras por pocillo. Los IDs de muestras se pueden mover y copiar.

Cuando se seleccione un método, en el lado izquierdo se mostrará una **Lista de todas las muestras desde diseño de placa**. En la lista solamente aparecen las muestras sin réplicas. Los IDs de muestras se pueden escribir en la lista o en la placa. Todos los cambios se muestran inmediatamente.

Botones de la barra de herramientas

Botón Ajustes	Este botón abre el cuadro de diálogo Configuración de listas de IDs de muestras , donde puede seleccionarse el número de IDs de muestras por pocillo.
Botón Editar	El botón Editar abre el cuadro de diálogo Editar pocillo . Se pueden editar los ID de muestras del pocillo marcado. Siga leyendo para obtener más información.
Botón Llenar automáticamente	Pueden seleccionarse la placa completa o una parte de la placa para completarse automáticamente con IDs de muestras. Después de seleccionar los pocillos correspondientes, haga clic en Llenar automáticamente para abrir el cuadro de diálogo Selección de llenado automático. Siga leyendo para obtener más información.
Botón Eliminar	Se puede seleccionar la placa completa o parte de la placa. Esta función borra los ID de muestras de los pocillos seleccionados inmediatamente. Ningún cuadro de diálogo pide confirmación. Haga clic en Borrar o pulse la tecla Supr para borrar los ID de muestras de los pocillos seleccionados.
Botón Deshacer	Deshace la última acción. Haga clic en Deshacer o pulse Ctrl-Z para anular una acción previa.

Botón Rehacer	Repite la acción que se ejecutó antes del paso de deshacer. Haga clic en Rehacer o pulse Ctrl+Y para repetir una acción que se había anulado.
Formato de impresión	Seleccione el diseño para imprimir la lista de IDs de muestras: • Tabla horizontal • Tabla vertical • Matriz
Presentación preliminar	Abre la presentación preliminar de la lista de IDs de muestras impresas.
Botón 10%	Fija la vista del diseño de placas al 10%.
Botón 100%	Fija la vista del diseño de placas al 100%.
Casilla de verificación Zoom	En caso de que la placa de pocillos conste de un gran número de pocillos, las asignaciones de pocillos individuales no se pueden mostrar en el modo de zoom al 100%. En ese caso, el botón Modo zoom permite un zoom en el área marcada. Cuando se activa, el usuario puede seleccionar un área para ampliarla haciendo clic y arrastrando un marco sobre el área de diseño deseada. Haga clic con el botón secundario del ratón para reducir hasta el 100%.

Menú contextual del Diseño de placa

Si se hace clic con el botón secundario en el diseño de placas, se mostrará un menú contextual para los pocillos marcados. Están disponibles los siguientes comandos:

Menú	Descripción		
Resumen	Se muestra el cuadro de diálogo Resumen.		
	Siga leyendo para obtener más información.		
Cortar	Los ID de muestras de los pocillos marcados se borran de los pocillos y se copian al portapapeles.		
Copiar	Los ID de muestras de los pocillos marcados se copian al portapapeles.		
Pegar	Los ID de muestras copiados al portapapeles se pegan en los pocillos marcados.		
Borrar	Haga clic en Borrar o pulse la tecla Supr para borrar los ID de muestras de los pocillos marcados.		
Copiar en formato ASCII	Los ID de muestras de los pocillos marcados se copian al portapapeles en formato ASCII. De este modo, se pueden pegar fácilmente en otras aplicaciones (p. ej., Excel, Bloc de notas).		
Pegar en formato ASCII	Los ID de muestras se pueden pegar del portapapeles en formato de archivo ASCII desde otras aplicaciones. Siga leyendo para obtener más información.		
Editar	La opción Editar abre el cuadro de diálogo Editar pocillo . Se pueden editar el ID de muestra del pocillo marcado. Siga leyendo para obtener más información.		



Menú	Descripción
Selección de Ilenado automático	Pueden seleccionarse la placa completa o una parte de la placa para completarse automáticamente con IDs de muestras. Después de seleccionar los pocillos correspondientes, al seleccionar la opción Llenar automáticamente se muestra el cuadro de diálogo Selección de llenado automático . Siga leyendo para obtener más información.



Aviso

Al editar una lista de IDs de muestras, puede arrastrar y colocar. Utilice el botón principal del ratón para seleccionar y arrastrar elementos.

Editar pocillo...

Se puede acceder al cuadro de diálogo **Editar pocillo** a través del menú contextual o el botón de la barra de herramientas.

Se pueden definir IDs de muestras para todos los pocillos de la microplaca.

Se puede utilizar un máximo de tres ID de muestras por pocillo.

Los ID de muestras deseados para el pocillo marcado se deben introducir en el cuadro de diálogo **Editar pocillo**. Esto se debe repetir para cada pocillo que se rellenará con IDs de muestras. Para trabajar con más comodidad se pueden utilizar las funciones de navegación del cuadro de diálogo **Editar pocillo**. Haga clic en **Aceptar** para cerrar el cuadro de diálogo. Se guardarán los ID de muestras para el pocillo marcado más reciente.

El cuadro de diálogo Editar pocillo contiene los siguientes elementos:

Botones Izquierda, Derecha, Arriba y abajo	Los botones de dirección se pueden usar para seleccionar otro pocillo en la placa mientras se deja el cuadro de diálogo abierto y se guardan los cambios del pocillo actual. Las Teclas de flecha también se pueden utilizar para este tipo de navegación.
Casilla de verificación Tras la entrada mover en sentido vertical	Al hacer clic en Intro o en Tabulador , el campo de entrada se mueve al siguiente pocillo. Use la casilla de verificación para especificar la dirección del movimiento, verticalmente u horizontalmente a través de la placa.
Campos de texto Editar ID de muestra	En cada pocillo se pueden escribir hasta tres IDs de muestras. Por lo tanto, la lista contiene tres campos de texto ID de muestra y una la casilla de verificación Autoincremento asociada a cada uno de ellos.
	En los campos de texto se pueden escribir los ID de muestras para el pocillo marcado. Los ID de muestras pueden contener letras, números o una combinación de ambos. Por ejemplo: Smp 1. Se pueden introducir hasta 100 caracteres alfanuméricos. Se puede rellenar toda la placa con IDs de muestras.



TECAN.

Selección de llenado automático

Se puede acceder al cuadro de diálogo **Selección de llenado automático** a través del menú contextual o el botón de la barra de herramientas.

Este cuadro de diálogo permite al usuario llenar automáticamente con IDs de muestras los pocillos seleccionados.

En el cuadro de diálogo **Selección de llenado automático** seleccione la placa completa o partes de la placa y escriba los IDs de muestras que desee y los ajustes de incrementos para los pocillos marcados. Haga clic en **Aceptar** para cerrar el cuadro de diálogo y llenar los pocillos seleccionados con los IDs de muestras definidos. En cada pocillo se pueden escribir hasta tres IDs de muestras. Por lo tanto, la lista contiene tres campos de texto **ID de muestra**, una casilla de verificación **Autoincremento**, un campo de texto **Índice inicial** y una casilla de verificación **Vertical** asociada a cada uno de ellos.

El cuadro de diálogo **Selección de llenado automático** contiene los siguientes elementos:

Casilla de verificación Utilizar ID	Solamente se puede tener acceso a la casilla de verificación Utilizar ID si se seleccionan números idénticos de IDs de muestras. La marca de comprobación determina si se aplicará la función de autoincremento al ID de muestra seleccionado.
Campos de texto ID de muestra	Los IDs de muestras para el pocillo marcado se puede escribir en los campos de texto ID de muestra . Los IDs de muestras pueden contener letras, números o una combinación de letras y números.
Casilla de verificación Autoincremento	Se debe seleccionar la casilla de verificación Autoincremento si se deben ser numerar automáticamente (incremento = 1) los IDs de muestra de los pocillos.
Campo de texto Índice inicial	El campo de texto Índice inicial solamente se puede editar si está seleccionada la casilla de verificación Autoincremento . Se debe escribir el valor de partida para el incremento automatizado de los ID de muestra.
Casilla de verificación Vertical	A la casilla de verificación Vertical solamente se puede acceder si está seleccionada la casilla Autoincremento . Si está seleccionada esa casilla, los IDs de muestras se escribirán y numerarán en dirección vertical. Si está desactivada, los IDs de muestras se escribirán y numerarán en dirección horizontal.



Resumen de pocillos

Se puede tener acceso cuadro de diálogo **Resumen de pocillos** mediante el menú contextual.

Este cuadro de diálogo muestra la perspectiva general de todos los IDs de muestras definidos y los ajustes relacionados con la lista de IDs de muestras de un pocillo seleccionado.

El cuadro de diálogo Resumen contiene los siguientes elementos:

Árbol de información	El árbol de información ofrece información general para el pocillo seleccionado sobre todos los parámetros definidos para pocillos. El árbol se rellena según la información disponible. Para las listas de IDs de muestras, se muestran los IDs de muestras, el estado de pipeteado, etc.
Botones Izquierda, Derecha, Arriba y Abajo	Los botones de dirección se pueden usar para seleccionar otro pocillo en la placa mientras se deja el cuadro de diálogo abierto.
Expandir todos / Reducir todos	Este botón muestra el árbol de información expandido hasta el nivel superior o contraído hasta el primer nivel.

Pegar en formato ASCII

Se puede tener acceso a la función **Pegar en formato ASCII** mediante el menú contextual.

Esta función permite transferir datos desde otras aplicaciones (por ejemplo: Excel, Bloc de notas) al diseño de placa de **magellan**.

En primer lugar hay que seleccionar en el otro programa los datos relevantes y copiarlos al portapapeles. Las filas deben estar separadas por avances de línea y las columnas por tabulaciones. Esto se hace automáticamente en Excel cuando se seleccionan varias celdas.

La función **Pegar en formato ASCII** pega los datos copiados en formato ASCII del portapapeles a los pocillos. Los datos siempre se insertan empezando por la posición A1.



5.2.3 Importar lista de IDs de muestras

Las listas de IDs de muestras creadas con otros programas o las listas de IDs de muestras definidas con otros formularios de equipos analíticos deben ser importadas a **magellan** para poder utilizarlas.



PRECAUCIÓN

ALGUNAS LISTAS DE IDS DE MUESTRAS PUEDEN CONTENER INFORMACIÓN ADICIONAL, COMO CONCENTRACIONES Y DILUCIONES, QUE PODRÍAN AFECTAR A LA EVALUACIÓN DE LOS DATOS DE MEDICIÓN.

Haga clic en el botón **Importar** del cuadro de diálogo **Configuración de listas de IDs de muestras** examinar los archivos externos de IDs de muestras existentes. Se mostrará el cuadro de diálogo **Abrir** de Windows[™].

Esa ventana ofrece una lista de **Nombres de archivo** que corresponden al **Tipo de archivo** seleccionado. El **Tipo de archivo** se puede seleccionar en una lista desplegable asociada.

Se admiten los siguientes tipos de archivos externos:

- · Archivos Easy .esy
- · Archivos Tecan .tpl
- Archivos DD1 .dd1
- Archivos Hamilton .pro
- Archivos APL .apl
- Archivos Gemini .csv
- Archivos de formato personalizado .txt

Seleccione el archivo solicitado para importarlo y haga clic en el botón **Abrir**. Se cerrará el cuadro de diálogo **Abrir** de Windows[™] y se cargará el archivo.

Tras confirmar el cuadro de diálogo **Configuración de listas de IDs de muestras**, con **Siguiente** se abrirá la ventana de placas **Importar o modificar lista de IDs de muestras** que contiene las IDs de muestras.

Al botón **Importar** también se puede acceder mediante el cuadro de diálogo **Insertar lista de IDs de muestras**, que está disponible en la ventana **Iniciar medición** (botón **Insertar**) y vista de placa **Evaluar resultados** (menú de barra de herramientas **Edición**).

STOP



ASEGÚRESE DE SELECCIONAR LA LISTA DE IDS CORRECTA, DE LO CONTRARIO LOS RESULTADOS SE ASIGNARÁN A OTRA LISTA.



Aviso Al importar listas de IDs de muestras en formato .tpl, .dd1 o personalizado, se puede incluir información adicional que modifica las definiciones de evaluación.



Aviso Cuando se importan listas de IDs de muestras con diluciones, la dilución no debe ser mayor que 1:1e+12 (magellan Standard y magellan Tracker).



Archivos Easy .esy

Estos archivos constan de líneas que contienen IDs de muestras y el nombre del pocillo.

nombre de pocillo:	alfanumérico, 3 caracteres máx.
ID de	alfanumérico, 20 caracteres máx.

muestra:

Las líneas están separadas por CR + LF.

Ejemplo

C01

B01

A01 0001

H02 0002

G02 0003

F02 0004

E02 0005

Interpretación

Cuando magellan procesa un archivo .esy

- El número de IDs de muestras se ajusta a 1
- El ID de muestra se establece como el ID de muestra en el pocillo

Archivos Tecan .tpl

Estos archivos (<ID_de_placa>.TPL) constan de líneas que contienen parámetros separados por punto y coma. El formato es:

H;fecha;hora

D;núm.test;ID.muestra + nombre.control;posición;predilución;errores

L;

El archivo empieza por H y termina en L. Los campos se definen como sigue:

fecha;hora;núm.test;ID.muest ra	igual que en la lista de trabajos (BES3)
nombre del control	alfanumérico, 8 caracteres máx. ¡Aviso! Esta entrada contiene el ID de muestra o el nombre de un control
posición	alfanumérico, 3 caracteres máx.
predilución	numérico, 8 caracteres máx. más 1 decimal (separado por un '.')

errores

alfanumérico, 3 caracteres máx. Si se produjo un error durante el pipeteado, aquí se muestra uno de los siguientes caracteres: Primer carácter		
<espacio> no hay error de líquido</espacio>		
L:	no hay líquido o no es suficiente	
Segundo carácter		
<espacio></espacio>	no hay error de código de barras	
M: manualmen	código de barras introducido te	
Tercer carácter		
<espacio></espacio>	no hay tiempo sobrepasado	
O:	tiempo sobrepasado	

Las líneas se separan con CR + LF.

Ejemplo

H;29-02-94;12:47:11

D;1;NEG;A1;1.0;

D;1;POS;B1;1.0;

D;1;0001;C1;1.0

L;

Interpretación

Cuando magellan procesa un archivo .tpl

- El número de IDs de muestras se ajusta a 1
- El ID de muestra o el nombre de control se establece como el ID de muestra en el pocillo
- Se guardan indicadores de error para cada pocillo
- · La predilución se utiliza como la dilución de la posición del pocillo

Archivos DD1 .dd1

Estos archivos constan de líneas que contienen parámetros de longitud fija.

El formato es:

Primera línea: Fecha (DD-MM-AA)

Segunda línea: Hora (HH:MM:SS)

Campo	Pos.	Longitud	Formato / Descripción
TestNr	1-3	3	Numérico (001-049)
TestName	4-23	20	Alfanumérico
SampleID	24-35	12	Alfanumérico
ControlName	36-43	8	Alfanumérico
PlateID	44-51	8	Alfanumérico
PlatePos	52-55	4	Alfanumérico
Errors	56-61	6	Alfanumérico

Las líneas se separan con CR + LF.



Ejemplo

13-04-1999			
11:01:32			
953test1 KBR		KBR3	H1
953test1 KBR		KBR3	G1
953test1 KBR		KBR3	C1
953test1 KBR		KBR3	B1
953test1 KBR		KBR3	A1
953test1 KBR	0001	KBR3	H2
953test1 KBR	0001	KBR3	G2
953test1 KBR	0001	KBR3	F2
953test1 KBR	0001	KBR3	E2
953test1 KBR	0001	KBR3	D2

Interpretación

Cuando magellan procesa un archivo .dd1

- El número de IDs de muestras se ajusta a 1
- · <Testname> especifica el nombre del método de magellan en el pocillo
- <PlatePos> especifica el pocillo
- · <SampleID> o <ControlName> especifica el ID de muestra de magellan
- <Errors> especifica el marcador de pipeteado: L: Error de líquido, C: Error de coágulo, E: Introducido manualmente

Cuando se lee un archivo .dd1 se modifica el diseño actual utilizando el diseño definido en los distintos métodos, donde los nombres se especifican en el archivo .dd1.

Archivos Hamilton .pro

Estos archivos constan de líneas que contienen el ID de muestra y el nombre del pocillo.

También hay disponible una información de encabezado.

Nombre de pocillo: alfanumérico, 3 caracteres máx.

ID de muestra: alfanumérico, 20 caracteres máx.

Las líneas se separan con CR + LF.

Ejemplo

14ESP-AK-Bommeli	-> Encabezado
08403009070	-> Encabezado
SendInfoHost=00	-> Encabezado
SendInfoPrnt=00	-> Encabezado
SendInfoDisk=00	-> Encabezado



A01091111111110

A02090325228840

A03090325228930

A04090325235730

A05090336975200

Interpretación

Cuando magellan procesa un archivo .pro

- El número de IDs de muestras se ajusta a 1
- El ID de muestra se establece como el ID de muestra en el pocillo
- · No se tiene en cuenta la información del encabezado

Archivos APL .apl

5.30 o posterio	r. El formato	o de archivo	se define como sigue:
Campo	Pos.	Longitud	Formato / Descripción
PlateID	1-8	8	Alfanumérico
PlatePos	9-12	4	Alfanumérico
TestNumber	13-15	3	Numérico
TestName	16-27	12	Alfanumérico
SampleID	28-39	12	Alfanumérico
ControlName	40-47	8	Alfanumérico
Result_1	48-55	8	Alfanumérico
Result_2	56-63	8	Alfanumérico
Errors	64-69	6	Alfanumérico
Date	70-79	10	Alfanumérico
Time	80-87	8	Alfanumérico
CutOff	88-95	8	Alfanumérico

Descripción de archivos .apl creados por MiniSwift 1.3 o posterior o Minilyser 5.30 o posterior. El formato de archivo se define como sigue:

Las líneas se separan con CR + LF.

Ejemplo

990520S1A1	2	test1_2	BLANK		05-20-199919:12:53
990520S1B1	2	test1_2	BLANK		05-20-199919:12:53
990520S1C1	2	test1_2	NC1	LR	05-20-199919:12:53
990520S1D1	2	test1_2	PC1		05-20-199919:12:53
990520S1E1	2	test1_2	0000000010		05-20-199919:12:53
990520S1F1	2	test1_2	0000000010	V	05-20-199919:12:53

Interpretación

Cuando magellan procesa un archivo .apl

- El número de IDs de muestras se ajusta a 3
- <Testname> especifica el nombre del método de magellan en el pocillo
- <PlatePos> especifica el pocillo
- <SampleID> especifica el ID de muestra 1 de magellan



- <ControlName> especifica el ID de muestra 2 de magellan
- <Errors> especifica el ID de muestra 3 de magellan
- <Errors> especifica el marcador de pipeteado: los errores de pipeteado L, B, N en la posición 1 no son críticos para las muestras, pero son críticos para los controles. La verificación de lectura y los errores de código de barras tampoco son críticos.

Archivos Gemini .csv

Descripción de archivos .csv creados por Gemini 3.40 y 3.60. A continuación se muestra un ejemplo de un archivo de salida en formato Excel:

Position	Tube ID	ScanError	SRCRack	SRCPos	SRCTubeID	Volume
Error	SRCRackID	GridPos	SiteOnGrid	TipNumber	DetectVol	Time
	Version 2.00	3091	Microplate	12	8	
0	9900001	18	1			3.8.1998
1		0	Tube 13*85mm	1	0001	10
0	013_000064	1	1	1		11:55:42
3		0	Tube 13*85mm	2	0002	20
0	013_000064	1	1	2		11:55:42
5		0	Tube 13*85mm	3	0003	30
0	013_000064	1	1	3		11:55:42
7		0	Tube 13*85mm	4	0004	40
0	013_000064	1	1	4		11:55:42

El formato de archivo se define como sigue:

La primera línea contiene los encabezados de columna, que son los mismos para todos los formatos, excepto para algunos que solo admiten caracteres en mayúsculas. La línea segunda muestra información acerca del rack o del rack de bandas para el cual fue creado el archivo. Solo son válidas las columnas Tube ID, SRCRack, SRCPos, SRCTubeID, Error, SRCRackID, GridPos, SiteOnGrid y Time.

Con Gemini 3.60, SRCPos contiene el número de pocillos en el eje x, SRCTubeID contiene el número de pocillos en el eje y Volume contiene la línea que se va a leer.

La tercera y las líneas que siguen son entradas para los pocillos utilizados de un rack o los tubos utilizados de un rack de bandas.

Las columnas contienen la siguiente información (Columna\Rango\Explicación):

- Posición\1... número de pocillos\Posición del pocillo en el rack. La secuencia de numeración es de atrás a adelante y después de izquierda a derecha.
 (1 = A1, 2 = B1, ...) Si la microplaca ha sido pipeteada en modo vertical (p. ej., dimensiones 8x12, iso. 12x8), la secuencia de numeración es de izquierda a derecha y después de adelante a atrás. (1 = H1, 2 = H2, ...)
- TubeID\Max. 32 caracteres\Esta columna solamente se rellena para racks de bandas y cuando se escanean los códigos de barras de los tubos con un comando POSID. En la fila 2, esta columna muestra la versión de Gemini.

 ScanError\0... 4294967295\Esta columna contiene posibles errores de escaneado. Tenga en cuenta la siguiente tabla con los códigos de error. En la fila 2, esta columna muestra el número de serie de Genesis.

TECAN.

- SRCRack\Max. 32 caracteres\Aquí Gemini notifica el nombre de configuración del rack fuente para los comandos de dispensar. En la fila 2, esta columna muestra el nombre de configuración del rack para el cual se escribió el archivo de salida.
- SRCPos\1 .. número de pocillos\Esta columna indica la posición en el rack de fuente de dónde llega el líquido para los comandos de dispensar.
- SRCTubeID\Max. 32 caracteres\Si la fuente para un comando de dispensar fuera un rack de bandas, Gemini aquí notifica el código de barras del tubo de origen.
- Volume\-7158278 .. +7158278\Esta columna contiene el volumen, en μl, que fue dispensado (valor positivo) o aspirado (valor negativo).
- Error\0 .. 4294967295\Esta columna contiene posibles errores de pipeteado. Tenga en cuenta la siguiente tabla con los códigos de error. En la fila 2, esta columna muestra los errores de escaneado.
- SRCRackID\max. 32 caracteres\Para los comandos de dispensar, este es el código de barras del rack de origen. En la fila 2, esta columna muestra el código de barras del rack, para el cual se escribió el archivo de salida.
- GridPos\1..67 Esta es la posición de retícula del rack de origen para los comandos de dispensar.
- SiteOnGrid\1 .. 128\Aquí Gemini muestra la posición y (posición en el portador) del rack de origen para los comandos de dispensar.
- TipNumber\1 .. 8\Aquí Gemini indica el número de punta utilizada para pipetear en/desde este pocillo.
- DetectVol\0 .. +7158278\Si se utiliza detección de líquidos, aquí se indica el volumen detectado en el pocillo antes del comando de dispensar o aspirar.
- Time\...\Esta es la hora (en formato hh:mm:ss) a la que se utilizó el pocillo.

Las líneas están separadas por CR + LF. Las columnas se separan por comas o por punto y coma. El separador de columnas se detecta según la primera línea de encabezado.

Interpretación

Se cuenta el número de líneas separadas por caracteres de nueva línea. Este número se reduce por 2 y se usa para definir el formato de placa, que puede ser 8x12, 16x24, 24x36, 32x48 (filas x columnas).

Con Gemini 3.60, la información de formato se extrae de la segunda línea, SRCPos y SRCTubeID. Se permite usar varias líneas por pocillo. El número de la línea segunda, Volume, define la línea que se va a utilizar. Si no existe la línea indicada, no se podrá leer la lista de IDs de muestras. Si alguna línea contiene un número de error mayor que cero, en función del número de error se establece como error crítico o error no crítico.

- Posición: Posición del pocillo (p. ej., 1 = A1, 2 = B1, ...)
- Tube ID: se ignora
- Scan error: si el número de error es mayor que cero, se establece un error crítico
- SRCRack: ID de muestra 2
- SRCPos: si SRCTubeID está vacío, se utiliza como ID de muestra 1 (SRCRackID + "_" + SRCPos)
- SRCTubeID: si no está vacío, se utiliza como ID de muestra 1
- Volume: se ignora



- Error: ID de muestra 3, si el número de error es mayor que cero, se establece un error crítico
- SRCRackID: si SRCTubeID está vacío, se utiliza como ID de muestra 1 (SRCRackID + "_" + SRCPos)
- GridPos: se ignora
- TipNumber: se ignora
- DetectVol: se ignora
- Time: se ignora

Archivos de formato especificado por el usuario *.*

Al importar archivos de formato personalizado se muestra el cuadro de diálogo Formato especificado por el usuario. El cuadro de diálogo Formato especificado por el usuario contiene los siguientes elementos:

Cuadro de grupo Tipo de placa	En el cuadro de grupo Tipo de placa se puede especificar el formato de la placa. Si el formato ya está especificado, se desactivan los campos de edición.
Filas	Introduzca el número de filas.
Columnas	Introduzca el número de columnas.
Cuadro de grupo Formato de archivo	En el Cuadro de grupo Formato de archivo se puede definir el formato del archivo especificado.
Extensión de archivo	En el campo Extensión de archivo se muestra la extensión del archivo y el campo está desactivado.



Aviso

Cuando se utiliza esta función en el asistente para Crear/Editar un método, el campo de extensión de archivo está habilitado y es preciso especificar la extensión de archivo.

Cuadro combinado Separador columna	Seleccione el separador de columnas utilizado en el archivo. El separador de fila debe ser CR+LF.
Empezar en la línea	Introduzca un número mayor que 1 si desea ignorar las líneas de encabezado.

Ejemplo

Usamos la siguiente configuración:

Tipo de placa

- 2 filas
- 3 columnas

Formato

- separador de columnas = Tabulador
- empezar en la línea = 2
- columna 1 = Posición de placa (A1, A2, ...)
- columna 2 = ID de muestra 1
- columna 3 = estado de pipeteado



Ejemplo:		
Archivo personalizado de Tecan		
A1	BLANK	0
A2	NC	0
A3	SM001	0
B1	SM002	3
B2	SM003	0
B3	SM004	0

5.2.4 Guardar listas de IDs de muestras

Haga clic en **Siguiente** en la ventana **Importar o modificar una lista de IDs de muestras** y se mostrará la ventana **Guardar como**, donde se puede guardar la lista creada o modificada de IDs de muestras:

Campo Nombre de archivo	Este campo sirve para escribir o modificar el nombre de archivo de la lista de IDs de muestras (.smp).
Campo Comentarios del archivo	El campo Comentarios del archivo permite añadir algunos comentarios al archivo que se mostrarán, por ejemplo, en los cuadros de diálogo Abrir archivo .
Campo Comentarios del registro de modificaciones	El campo Comentarios del registro de modificaciones permite añadir algunos comentarios, que se guardarán en el registro de modificaciones (solamente disponible para magellan Tracker).
Botón Registro de modificaciones	Muestra el registro de modificaciones de la lista de IDs de muestras (solamente disponible en magellan Tracker). Para obtener información adicional, consulte 4.4 Guardar métodos - Registro de modificaciones.

El botón **Guardar y finalizar** solamente estará habilitado cuando se haya modificado un documento. Ese documento modificado se puede guardar con su nombre de archivo actual o con un nombre diferente.

Haga clic en **Guardar y finalizar** para guardar la lista de IDs de muestras y cerrar el asistente **Crear o editar lista de IDs de muestras**.



6. Asistente Iniciar medición

6.1 Introducción

Para abrir este asistente, en la ventana Lista de asistentes, haga clic en Iniciar

medición. Haga clic en Continuar en la página de bienvenida y se mostrará el cuadro de diálogo Seleccionar un archivo:

Seleccionar un archivo						\times
O Obtener datos sin procesar						
 Usar método previamente definido 			Nostrar:	Archivos de este i	nstrumento 🗸	
O Activar favorito						
nth		Name	Notas		Estado	
Anuda						
			HAGA	SU SELECCIÓ	N 🕨	
Cancelar <<< Atras						
				© 2021 Tecan		
Obtener datos sin procesar	Se utiliz mediant	a para generar datos e re el ajuste de los parár	n bruto metros	rápidamer de medició	nte	
	necesar	ios.				
Usar método predefinido	Se emp ya defin	lea para realizar medic idos.	iones b	basadas en	métodos	
Activar favorito	Se usa uno de l	para seleccionar en la l los métodos usados co	lista de n más	íconos en frecuencia.	umerados	5

6.2 Obtener datos sin procesar

6.2.1 Obtener datos sin procesar con el instrumento SUNRISE

Seleccione **Obtener datos sin procesar** y haga clic en **Siguiente** para abrir el cuadro de diálogo **Parámetros de medición**.

En el cuadro de diálogo **Parámetros de medición** se ofrecen las siguientes opciones:

Medición de punto final	Utilice este botón para realizar una medición individual.
Medición cinética	Este tipo de medición se utiliza para mediciones múltiples dentro de un intervalo de tiempo especificado.
Parámetros de medición	Haga clic en este botón para definir los parámetros del cuadro de diálogo Parámetros de medición .



ADVERTENCIA

SE RECOMIENDA COMPROBAR TODOS LOS PARÁMETROS DE MEDICIÓN (INCLUYENDO AQUELLOS QUE NO SON EMPLEADOS EN LA MEDICIÓN ACTUAL) YA QUE LOS PARÁMETROS DEL MÉTODO DE MEDICIÓN PREVIO SEGUIRÁN TODAVÍA EN EFECTO.

6.2.2 Obtener datos sin procesar con el instrumento INFINITE F50

Cuando haya conectado un instrumento INFINITE F50, se mostrará la siguiente ventana después de seleccionar **Obtener datos sin procesar**:

Parametrus de medición - Urbeita F90			
Herramientas de laboratorio 💈	state and the second se		Selection
+ Rec	Definición de placa (GRESOR) - Greiner 94 Mail Transport	U Detailes	nacia pereccorsezo
Siccer of pace		Utilizar una parte de la placa	
T Heliciones 2	👆 🖛 Sección de placa	2	
4 Absorbencia			
Access 2	A 000000000000000000000000000000000000		
X Aptación			
🔨 Noverplace			
Cevitaca 2	r 0000000000		
G Cide enético	* 6990666669000		
Se Condición covitica			
Otras opciones #	4 - Absorbencia	3	
🔁 Conetana	Longitud de onda Etiqueta		
 Solicitud del usania Esperar (trespontration) 	Hediośni 405 twi v Norbini: jubel v		
😸 äncubación	R relative interaction and the second		
Número de placas: [32]			
Ayuda		Activate Wind	lows
Cancelar <<< Atrās		NELECCIONAR PARAMETROS DE	MEDICION COM

Defina los parámetros de medición según la lista del lado izquierdo.

Para terminar la definición, haga clic en Seleccionar parámetros de medición

Para obtener información adicional, consulte las Instrucciones de uso del i-control.



6.3 Usar método predefinido

Haga clic en **Usar método predefinido** y seleccione un método en la lista **Nombre de archivo**.

Los métodos disponibles se muestran según las opciones de la lista desplegable Mostrar, es decir, Todos los archivos, Archivos de este instrumento, Mis archivos, Archivos firmados o Últimos métodos seleccionados.

6.4 Activar favorito

Haga clic en **Activar favorito** para seleccionar en la lista de iconos numerados un método predefinido como favorito (un método que se utiliza con frecuencia).

Haga clic en **Organizar favoritos** para añadir o quitar métodos de la lista de favoritos.

Organizar favoritos

Los archivos de métodos se pueden arrastrar desde la lista **Método** hasta uno de los 15 íconos de la lista **Favoritos** para guardarlos como favoritos.

)rganizar favoritos					
Método:					
Nombre de archivo		Notas		Estado	
ELISA_Qualitative_M1000.mth				No firmado	
Distance of the second	M1000.mth			No firmado	Añadir Eliminar
Métodos favoritos:	2	3	4	5	<u>^</u>
<vacío></vacío>	<vacío></vacío>	<vacío></vacío>	<vacío></vacío>	<vacío></vacío>	
46	47	8	49	40	
<vacío></vacío>	<vacío></vacío>	<vacío></vacío>	<vacío></vacío>	<vacío></vacío>	~
Instrumento favorito infi ELISA_Qualitative_M10	nite M1000 100.mth	Seleccior	nar	Aceptar	Ayuda Cancelar

El cuadro de diálogo Organizar favoritos consta de los siguientes elementos:

Lista Métodos	Enumera todos los métodos disponibles. Muestra el nombre del archivo y, si se han escrito previamente, las anotaciones correspondientes.
Botón ñ adir	Después de seleccionar un método en la lista Nombres de archivos de métodos y seleccionar un icono de método, haga clic en Añadir para añadir el método seleccionado a Favoritos .
Botón Eliminar	Seleccione un icono de método y haga clic en Eliminar para eliminar un método de Favoritos .

6.5 Iniciar medición con un método predefinido o favorito

Haga clic en **Haga su selección** respectivamente **Mostrar resumen** en el cuadro de diálogo **Seleccionar un archivo** para abrir el cuadro de diálogo **Iniciar medición**. Antes de que el instrumento empiece a medir, el usuario tiene que (si lo ha definido) contestar a las preguntas del informe o confirmar/cambiar las constantes.

Cuadro de diálogo Editar indicaciones

Si en el método se especificaron indicaciones al usuario, se mostrará el cuadro de diálogo **Editar indicaciones al usuario**:

Lista Respuesta a indicación	En cada campo de texto de esta lista se puede escribir cualquier texto de indicación al usuario.
Casilla de verificación Req.	La casilla de verificación Requerida junto a cada campo de texto Respuesta a indicación determina si la medición solamente podrá empezar si se ha escrito un texto.

Haga clic en **Continuar** para cerrar el cuadro de diálogo de indicaciones al usuario; se mostrará la ventana **Iniciar medición**. Si así se estableció en el método, se mostrará el cuadro de diálogo **Editar constantes**.

Cuadro de diálogo Editar constantes

El cuadro de diálogo **Editar constantes** solamente se mostrará si está activado el marcador Requerida para al menos una constante en el método seleccionado.

Lista Constantes	En cada campo de texto de esta lista se puede modificar el valor de la constante o el comentario.
Casilla de verificación Req.	La casilla de verificación Requerido junto a un campo de texto Valor constante indica que ese valor constante se debe confirmar. Para ello, siga las instrucciones de la página siguiente.

Si se han modificado las constantes de un método, los nuevos valores se pueden transferir y guardar automáticamente en el método correspondiente. Haga clic en **Modificar constantes de Método** para modificar el método actual y aplicar las nuevas constantes definidas. Tenga en cuenta que ese cuadro de diálogo solamente estará disponible para los usuarios de Magellan que tengan el derecho de usuario **Editar métodos**.

Haga clic en **Continuar** para cerrar el cuadro de diálogo de las constantes; se mostrará la ventana **Iniciar medición**.


i

Aviso

Si se utiliza magellan Tracker o magellan Standard con Administración de usuarios, un Operador de magellan predeterminado (consulte los capítulos 9.6 Derechos del usuario y 9.5.3 Cambiar usuario) solamente puede editar las constantes requeridas. Las constantes que no tengan activado el marcador Requerida aparecerán en gris y no se podrán editar.

Un Administrador predeterminado de magellan también puede editar las constantes sin el marcador Requerida activado.

En el caso de un Especialista en aplicaciones predeterminado de magellan depende del método que ejecute. Si utiliza un método no firmado, puede editar constantes con y sin el marcador Requerida. Si utiliza un método firmado, solamente puede editar las constantes con el marcador Requerida

Después de estas dos páginas opcionales, se muestra el cuadro de diálogo **Iniciar medición**:

Iniciar medición				×
Medición			Parámetros de medición	
Zona de trabajo:	07032018-008.wsp	Ciclo cin. arb	Modo de medición: Absorbencia	
Método:	Application Example_es.mt	Modificar diseño	Longitud de onda de referencia: 620 nm	
Lista de IDs de muestra:		Insertar	Modo lectura: normal	
Instrumento	/		Unidad: UD	
Utilizar apilador		Actual: 32.1 °C		
Introducir placa	Control de temperatura	Objetivo: 37.0 °C		
Movimientos	Control inyector 0	ptimiza la posición Z		
Tenga en cuenta lo sigui	ente:			
Aurida	/			
nyuud			INICIAR	
Cancelar	<< Atrás			

Cuadro de grupo Medición	Campo de texto Zona de trabajo : muestra el nombre predeterminado del archivo de zona de trabajo según se definió en Procesamiento automatizado de datos . Se puede cambiar el nombre de archivo de esta zona de trabajo.
	Botón Ciclo cin. arb : este botón está disponible si se realiza una medición cinética. Haga clic en este botón para mostrar el cuadro de diálogo Cinética de ciclo arbitraria. En este cuadro de diálogo las mediciones se pueden descomponer en diversas submediciones con distintos números de intervalos de tiempo y ciclos de medición. Cada una de esas submediciones se puede iniciar en cualquier momento, como por ejemplo: se puede guardar la zona de trabajo y continuar la medición otro día.

i



Esta opcio	ón no está disponible en instrumentos INFINITE F50.
	Campo de texto Método : se muestra el nombre de archivo del método seleccionado previamente.
	Botón Modificar diseño : en función de los derechos de usuario, permite modificar el diseño de las placas. Haga clic en el botón Modificar diseño para abrir el cuadro de diálogo Modificar diseño . El usuario podrá mover controles y retirar muestras, cambiar la concentración o modificar los parámetros de medición. Los cambios solamente se guardan en la zona de trabajo y no en el método seleccionado previamente. Este botón no estará disponible cuando se haya seleccionado Obtener datos sin procesar y después de insertar una lista de IDs de muestras.
	Campo de texto Lista de IDs de muestras : muestra el nombre de archivo de la lista de IDs de muestras cargada actualmente.
	Botón Insertar : consulte también las siguientes explicaciones. Haga clic en este botón para abrir el cuadro de diálogo Insertar lista de IDs de muestras . Después de seleccionar un archivo de lista de IDs de muestras, se muestra el cuadro de diálogo Lista de IDs de muestras con vista previa del diseño de placa para que el usuario confirme la selección.
Cuadro de grupo Instrumento	Las zonas de trabajo se guardarán con los nombres predeterminados de las zonas de trabajo. Después de que se hayan efectuado todas las mediciones de las placas, solo quedará abierta la primera zona de trabajo. Para abrir las restantes zonas de trabajo, utilice Abrir archivo o el asistente Evaluar resultados . No se puede utilizar el apilador mientras se realizan mediciones cinéticas .
	Botón Introducir o sacar placa: Haga clic para mover el portaplacas. El texto de este botón variará según la posición del portaplacas.
	Botón Movimientos : permite invocar movimientos adicionales del instrumento (placa de filtro, portaespejos, etc.)
	Campo de texto Temperatura actual : muestra la temperatura actual del instrumento.
	Campo de texto Temperatura objetivo : se muestra la temperatura objetivo que se define en el cuadro de diálogo Control de temperatura .
	Botón Control de temperatura : Haga clic en este botón para abrir el cuadro de diálogo Control de temperatura . Consulte el capítulo 3.1.2 Control de temperatura
Cuadro de grupo Parámetros de medición	Permite comprobar en un campo de texto todos los parámetros de medición configurados.



Cuadro de grupo Muestra información importante. Lea esa información y realice las acciones requeridas antes de aplicar el método. **Io siguiente**

Se mostrará una advertencia si el diseño está definido para pocillos que no se medirán.

Para empezar la medición, haga clic en Iniciar.

Si ejecuta un método multiplaca, se mostrará el cuadro de diálogo **Ejecutar medición** después de hacer clic en **Iniciar**.

jecutar medic	ón
Medición de	placa 🚺 de 3
Lista de IDs	de muestra:
-	
	Parar PAUSA EJECUTAR
	ntroducir placa

Contiene los siguientes elementos:

Medición de placa	Muestra el número de la placa procesada actualmente.
Lista de IDs de muestras	Inserte la lista de IDs de muestras apropiada para la placa seleccionada. Las listas de IDs de muestras se deben insertar placa a placa.



Aviso Solamente se pueden insertar listas de IDs de muestras creadas en magellan. No se admite la importación de otros formatos.

EJECUTAR	Haga clic en este botón para iniciar la medición.
PAUSA	Al hacer clic en Pausa se invocará el asistente Evaluar resultados; a continuación, se pueden anexar placas adicionales.
PARAR	Al hacer clic en Parar se invocará el asistente Evaluar resultados; no se pueden anexar placas a continuación.

Tras completarse la medición, aparecerá el cuadro de diálogo **Resultados** donde se muestran los resultados y los cálculos. Para obtener más información, consulte el capítulo 7 Asistente Evaluar resultados.



Insertar lista de IDs de muestras

La opción **Insertar lista de IDs de muestras** se utiliza para cargar en una zona de trabajo un archivo de lista de IDs de muestras definido previamente:

Campo Nombre de archivo	Asa lista muestra todos los archivos de listas de IDs de muestras disponibles (.smp). Haga clic en un nombre de archivo para seleccionar una lista de IDs de muestras.
Campo Notas	Mostrará los comentarios y las notas disponibles.
Botón Importar	Haga clic para abrir un cuadro de diálogo estándar Abrir de Windows. Permite importar archivos externos. Para obtener información adicional, consulte el capítulo 5.2.3 Importar lista de IDs de muestras.

Seleccione en el campo **Nombre de archivo** la lista de IDs de muestras que desee y haga clic en **Aceptar** para ver la lista de IDs de muestras en la ventana **Lista de IDs de muestras con vista previa del diseño de placa**. Este cuadro de diálogo muestra la lista de IDs de muestras junto con el diseño de placa definido. Si el método y la lista de IDs de muestras tienen formatos de placa diferentes, se mostrará un mensaje de error.

Cuando se selecciona la lista de IDs de muestras en la configuración de informe impreso, aparecerá en la copia impresa.

El cuadro de diálogo **Lista de IDs de muestras con vista previa del diseño de placa** muestra la lista de IDs de muestras junto con el diseño definido de la placa. Esto ayuda al usuario a asegurarse de que ha seleccionado la lista de IDs de muestras correcta. El cuadro de diálogo contiene los siguientes elementos:

Botón 10%	Haga clic para redimensionar el diseño de placa y mostrar al 10% el diseño de placa a pantalla completa.
Botón 100%	Haga clic para redimensionar el diseño de placa y mostrar al 100% el diseño de placa a pantalla completa.
Casilla de verificación Zoom	Cuando se activa esta función de Zoom , se pueden marcar los pocillos necesarios en el diseño de placa y los pocillos se ampliarán automáticamente para ajustarse a la pantalla completa.
Casilla de verificación Retire las muestras sin usar	Si se carga una lista de IDs de muestras solo rellenada en parte, el usuario puede seleccionar esta casilla de verificación para borrar del diseño todas las muestras no usadas. Si se ha definido dentro del diseño un control final (p. ej. en el pocillo H12), ese control se moverá automáticamente al pocillo después del último pocillo de muestra utilizado. De modo predeterminado, la opción Retire las muestras sin usar está seleccionada.

6.6 Estado de la medición

El cuadro de diálogo **Estado de la medición** se muestra durante la ejecución de la medición del instrumento y contiene los siguientes elementos:

Campo de texto Estado	Este campo de texto se actualiza durante la medición y contiene información de estado de la medición.
Cuadro de grupo Mostrar en pocillos	Seleccione el modo de representación de los datos en los pocillos: Valores : el valor medido se mostrará en el pocillo. Si se mide más de un valor se muestran todos los valores en un pocillo.
	Gráfico : esta opción estará habilitada si se puede mostrar un gráfico cinético. El gráfico se mostrará tan pronto como haya disponible más de un valor.
Botón Pausa	Para hacer una pausa en una medición cinética ejecutada en el siguiente momento posible, haga clic en el botón Pausa . En el cuadro de diálogo Haciendo una pausa en cinética haga clic en el botón Extraer placa / Introducir placa para mover dentro o fuera del instrumento el transporte de placa. Haga clic en el botón Continuar para proseguir con la ejecución cinética.
Botón Interrumpir	Haga clic en este botón para interrumpir la medición en el siguiente punto de tiempo disponible.
Pantalla Vista de placa	La pantalla Vista de placa muestra una visualización esquemática de la microplaca. Combinada con la opción Mostrar en pocillos permite personalizar el contenido de los pocillos. Los pocillos se actualizan tan pronto como el instrumento devuelve datos. Cuando se realiza la inyección o la distribución, el pocillo utilizado en ese momento se resalta automáticamente con
Campo de texto Ciclo	un color diferente. Si se realizada una medición cinética, el campo de texto Ciclo muestra el número del ciclo actual y el número máximo de ciclos cinéticos.
Pantalla Gráfico	Si se realiza una cinética o una medición de etiquetas múltiples, se mostrará en un tamaño más grande el gráfico cinético o de etiquetas múltiples del pocillo seleccionado actualmente a la derecha del cuadro de diálogo Estado de la medición .
Cuadro de grupo Ciclo previo	Mientras no comience la transmisión de los datos del ciclo actual, los campos de texto Datos sin procesar y Temperatura contienen, respectivamente, el valor y la temperatura del ciclo de medición anterior del pocillo seleccionado actualmente.
Cuadro de grupo Tiempo	El campo de texto Tiempo transcurrido muestra el período de tiempo desde el comienzo de la medición.
	El campo de texto Tiempo de ejecución estimado muestra el período de tiempo previsto para la medición completa.



Se puede cambiar la escala del eje de ordenadas del gráfico seleccionando o bien **Autoseleccionar rango** o **Seleccionar rango** (MÍN/MÁX).



Aviso

Se pueden seleccionar múltiples pocillos en la placa para ver las curvas cinéticas seleccionadas en un gráfico. Cuando se realizan mediciones cinéticas, los datos obtenidos se guardan automáticamente cada 30 minutos en el directorio de archivos de registro de magellan (consulte el capítulo 3.3 Archivos de registro).



7. Asistente Evaluar resultados

7.1 Introducción

El **asistente Evaluar resultados** está diseñado para ayudar al usuario a analizar los resultados de las mediciones. Permite visualizar los datos sin procesar, los datos de evaluación y los parámetros de evaluación, así como volver a evaluar los datos.

Resumen del flujo de trabajo

Inicie el **asistente Evaluar resultados** y seleccione una zona de trabajo existente para la evaluación (cuadro de diálogo **Seleccionar un archivo**)

Haga clic en **Haga su selección** para abrir el cuadro de diálogo **Resultados**, donde puede visualizar o imprimir los resultados (pestaña **Evaluar resultados**) y, si fuera necesario, modificar el método utilizado (pestaña **Editar método**).

Haga clic en **Finalizar** y utilice el cuadro de diálogo **Guardar** para guardar un archivo de zona de trabajo que contiene definiciones de métodos, datos del instrumento y listas de IDs de muestras (si se han especificado).

Los datos evaluados no se guardan en la zona de trabajo, sino que se calculados de nuevo cada vez que se abre la zona de trabajo.

7.2 Seleccionar un archivo

El cuadro de diálogo **Seleccionar un archivo** contiene los siguientes elementos:

Lista Nombre de archivo	La lista Nombre de archivo contiene la lista de todos los archivos (y las notas) dentro del directorio de la zona de trabajo estándar.
Cuadro combinado Mostrar	 En el cuadro combinado Mostrar se puede modificar la lista de archivos mostrados conforme a la selección: Todos los archivos Archivos de este instrumento Mis archivos: esta opción solo estará disponible si está habilitada la administración de usuarios (siempre está habilitada en magellan Tracker). Archivos firmados: solo disponible para magellan Tracker. Haga clic en el botón para crear definiciones personalizadas; consulte Criterios de selección de archivos a continuación. Archivos de ejemplo: solo estarán disponibles si se han instalado.

7.2.1 Criterios de selección de archivos

Todos los archivos Archivos de este instrumento	Aceptar Cancelar
Mis archivos	
O Archivos firmados	Ayuda
Archivo de zona de trabajo creado o r	nodificado
💽 Sin límite de tiempo	
ОНоу	
0	
🔘 La semana pasada	
C La semana pasada El mes pasado	

Haga clic en el botón para abrir el cuadro de diálogo **Criterios de selección de archivos**.

Seleccione una de las opciones bajo **Criterios de zona de trabajo** o escriba algunos caracteres en el campo de texto **Contiene** que se muestra en **Nombre de archivo**. Se puede seleccionar una de las opciones bajo **Archivo de zona de trabajo creado o modificado** para aplicar un filtro de tiempo o se puede escribir un número en el campo **Últimos ... días**. Haga clic en **Aceptar** para guardar esos ajustes.

Seleccione en la lista **Nombre de archivo** el archivo de zona de trabajo que desea evaluar y haga clic en **Haga su selección**.



Aviso

Si tiene un instrumento SUNRISE conectado pero selecciona un método de INFINITE F50, se mostrará el cuadro de diálogo Incongruencia de instrumento:

Conecte el instrumento adecuado

7.3 Ventana Vista de conjunto de la zona de trabajo

La ventana **Vista de conjunto de la zona de trabajo** muestra todos los datos disponibles de la zona de trabajo seleccionada actualmente.

Si el usuario que está trabajando actualmente tiene los derechos de magellan apropiados, también se puede hace clic en la pestaña **Editar método** para hacer algunos cambios en el método y volver a calcular los datos medidos.



La Vista de conjunto de la zona de trabajo está formada por los siguientes elementos:

- Ventana Diseño de placa: muestra los datos seleccionados para cada pocillo. Si se abre una zona de trabajo multiplaca, se mostrará en la esquina superior derecha un "contador" para cambiar entre las distintas placas.
- Barra de herramientas: para editar, hacer zoom, cambiar entre ciclos cinéticos y ver el resumen de un pocillo seleccionado.
- Ventana Información de datos: en esta ventana de cuadro de texto se muestran datos que no se pueden asignar a un pocillo individual (por ejemplo, parámetros de medición, definición de corte, estadística resultado, etc.).
- Pestaña Barra de control de Evaluar resultados: muestra todos los nombres de conjuntos de datos disponibles. La selección de un conjunto de datos mostrará los datos en el diseño de placa o mostrará los datos en la ventana Información de datos.
- Pestaña Barra de control de Editar método: muestra la configuración y los parámetros del método cargado.

El tamaño de estas tres últimas ventanas se puede cambiar por medio de las barra divisoras que separan las ventanas.



Diseño de placa

La vista muestra un diseño esquemático de una microplaca. Las filas se marcan alfabéticamente y las columnas numéricamente.

En la barra de control de la izquierda se muestran todos los datos disponibles. Los datos seleccionados aparecen en el diseño de placa, en el tablero con texto o en una ventana de gráfico aparte en función de los datos seleccionados.

Menú contextual de Diseño de placa

Para mostrar un resumen de todos los datos de un pocillo, seleccione el pocillo y seleccione **Resumen** en el menú del botón secundario del ratón.

Botones de la barra de herramientas

En los siguientes párrafos se describen los botones disponibles en la barra de herramientas.

En función de los datos seleccionados se muestran distintos botones: cuando se selecciona **Datos del instrumento** está disponible el botón **Editar**; cuando se selecciona **Concentración** o **Datos cinéticos**, está disponible el botón **Curva/Gráfico**.

Los siguientes botones solo están disponibles en el asistente Eval	uar
resultados.	

Editar	Haga clic en el botón Editar para editar u ocultar datos sin procesar.
Botón 🛄	Haga clic en este botón para ver el cuadro de diálogo Resumen .
Botón 10%	Haga clic en este botón para redimensionar el diseño de placa y mostrar el 10% del diseño de placa a pantalla completa.
Botón 100%	Haga clic en este botón para redimensionar el diseño de placa y mostrar el 100% del diseño de placa a pantalla completa.
Casilla de verificación Zoom	Cuando se selecciona la casilla de verificación Zoom se pueden marcar los pocillos necesarios en el diseño de placa y los pocillos etiquetados se ampliarán automáticamente para ajustarse a la pantalla completa.
Lista desplegable Ciclo de medición	Si la medición es una medición cinética, en la lista desplegable se puede seleccionar el ciclo de medición deseado.

Ventana de información de datos

Haga clic, por ejemplo, en datos de **Otras opciones** en la pestaña **Evaluar resultados** para ver la información correspondiente en la ventana **Información de datos**.

7.4 Pestaña Evaluar resultados

Todos los nombres de conjuntos de datos disponibles se muestran en la barra de control **Datos disponibles**. Hay varias opciones de visualización:

- Se muestran en la ventana de diseño de placa datos seleccionados que presentan un valor para cada pocillo. Se pueden ver como valores o como pseudocolores junto con una escala de colores.
- Los datos cinéticos se pueden ver como un pequeño gráfico en cada pocillo.
- Los datos seleccionados que presentan una colección de datos que no se puede asignar a un pocillo individual se muestran en la ventana de información de datos (p. ej., Criterios de validación QC).
- Si se selecciona un gráfico (curva estándar, cinética, cinética de enzima, series de dilución, etiquetas múltiples, espectros), se mostrará el cuadro de diálogo **Gráfico**.

Si se ha ejecutado una secuencia de medición que consta de mediciones independientes individuales, se mostrarán los resultados de cada medición al seleccionar en una lista desplegable de la barra de herramientas el número de medición correspondiente (p. ej., ciclos cinéticos en una ejecución cinética).

Si fuera necesario, se puede hacer clic en el botón **Editar** para modificar los resultados de la medición. Los datos manipulados se marcan con un signo ~.



ADVERTENCIA

ESTA FUNCIÓN SOLAMENTE ES PARA USO EN INVESTIGACIÓN, NUNCA SE DEBE UTILIZAR PARA PRUEBAS DIAGNÓSTICAS.

En cada pocillo del diseño de placa, se pueden mostrar tres líneas de datos. En función de los criterios seleccionados en el panel de control de la ventana izquierda, se muestran distintos valores.

7.4.1 Menús

Archivo	Exportar a archivo ASCII
	Exportar a archivo ASTM (LIS)
	Exportar rastreo de muestras
	Exportar a Excel
	Método Exportar
	Imprimir
	Presentación preliminar
	Configurar impresora
	Fuente de impresión
	QC de placa a placa
Edición	Copiar a Excel
	Copiar a formato ASCII
	Pegar de formato ASCII
	Insertar lista de IDs de muestras
	Volver a calcular con otro método
Instrumento	Movimientos
	Control de temperatura
	Iniciar medición



7.4.2 Menú de Barra de herramientas: Archivo

Exportar a archivo ASCII

Haga clic en **Exportar a archivo ASCII** para exportar los datos como archivos ASCII (.asc).

Antes de ejecutar el procedimiento de exportación, se deben seleccionar los datos en el cuadro de diálogo **Exportación de datos** mediante la pestaña **Editar método** \rightarrow **Procesamiento de datos** \rightarrow **Exportación de datos**.

Seleccione la carpeta donde desea guardar los datos exportados. Escriba un nombre de archivo (.asc).

Al hacer clic en **Guardar** en el cuadro de diálogo **Guardar como**, los conjuntos de datos especificados en el cuadro de diálogo **Exportación de datos** se guardarán como un archivo ASCII.

Exportar a archivo ASTM (LIS)

En el menú **Archivo**, haga clic en **Exportar a archivo ASTM** para exporta datos en formato ASTM compatible con LIS (Laboratory Information Systems).

Seleccione en el cuadro de diálogo **Exportación de datos** los datos que desea exportar a través de la pestaña **Editar método** \rightarrow **Procesamiento de datos** \rightarrow **Exportación de datos**.

Exportar rastreo de muestras

En el menú Archivo, haga clic en Exportar rastreo de muestras para exportar datos al sistema de rastreo de muestras. Seleccione en el cuadro de diálogo Exportación de datos los datos que desea exportar. Use la pestaña Editar método \rightarrow Procesamiento de datos \rightarrow Exportación de datos.

Exportar a Excel

Para exportar de datos a Excel, haga clic en Exportar a Excel.

Seleccione en el cuadro de diálogo **Exportación de datos** los datos que desea exportar a través de la **pestaña Editar método** \rightarrow **Procesamiento de datos** \rightarrow **Exportación de datos**.

Método Exportar

Haga clic en Método Exportar para exportar un método.

Seleccione la carpeta donde desea guardar el método y escriba el nombre de archivo que desee. Además, puede añadir algunos comentarios en el campo de texto **Notas**.

Imprimir

En el menú **Archivo**, haga clic en **Imprimir...** o pulse **MAYÚS-P** para imprimir los datos como haya definido en la configuración de **Informe impreso** desde el método. Se abrirá el cuadro de diálogo **Imprimir** estándar de Windows donde se puede seleccionar la impresora, el rango de páginas y el número de copias.

Presentación preliminar

Haga clic en **Presentación preliminar...** para ver exactamente el aspecto de la copia impresa sin necesidad de imprimir realmente el documento.

Configurar impresora

Haga clic en **Configurar impresora...** para cambiar la impresora actual o modificar la configuración de la impresora.

Se mostrará el cuadro de diálogo estándar **Configurar impresora** de Windows. Seleccione la impresora, el tamaño del papel y la orientación de la copia impresa.



Fuente de impresión

Haga clic en **Fuente de impresión...** para seleccionar los detalles de fuentes, estilos de fuente, tamaños de fuente, color de fuente, etc. para futuras copias impresas.

La configuración de estos ajustes no tendrá ningún efecto en la presentación en pantalla. Es recomendable no utilizar fuentes grandes para evitar errores y deformar la presentación de los datos.

QC de placa a placa

Cuando se configura el **QC de placa a placa** en la definición del método, **magellan** compara el valor medio actual observado de un control con la desviación media y estándar previamente definida para este Método. (Consulte el capítulo 4 Asistente Crear o editar método).

Gráfico Levy-Jennings

Haga clic en **QC de placa a placa** y se mostrará la ventana **Gráfico Levy-Jennings**. El gráfico Levy-Jennings muestra la tendencia de los valores de control.



Sobre los ejes de ordenadas, los valores de datos (datos sin procesar, concentraciones, etc.), tal como se establecen en la definición del método, se aplican sobre la escala de tiempo en días (ejes de abscisas).

Se muestran los valores de control junto al valor medio y las líneas de distancia de las desviaciones estándar (1s, 2s, 3s).

Cuando el cursor se coloca sobre un punto de medición, una información sobre herramientas (tooltip) muestra información adicional (nombre de la zona de trabajo, etc.). Utilice las casillas de verificación de los elementos individuales de la lista **Mostrar** para seleccionar los distintos controles que desea visualizar.

Intervalo de tiempo

Haga clic en **Intervalo de tiempo** para evaluar zonas de trabajo dentro de un intervalo de tiempo definido.



Seleccionar zonas de trabajo

Haga clic en **Habilitar/Deshabilitar** para activar o desactivar la evaluación de algunas zonas de trabajo.

Activar/Desactivar zonas de trabajo	×
21072009-004 MISD	
✓21072009-005.wsp	
✓21072009-006.wsp	
✓21072009-007.wsp	
✓21072009-008.wsp	
21072009-009.wsp	
Aceptar Cancelar Ayuda	

Hoja de datos

Haga clic en Ver resumen para abrir el cuadro de diálogo Hoja de datos:

Columna Zona de trabajo	Muestra la lista de los archivos de zona de trabajo evaluados. A la izquierda del nombre de la zona de trabajo, una marca de comprobación verde indica el cumplimiento y una cruz roja indica el fallo de la evaluación de la regla de Westgard [®] (para obtener información adicional acerca de las Reglas de Westgard , consulte las Instrucciones de uso de magellan).
Columna Control	Se enumeran los distintos controles.
Columna Valor	Se muestra el valor medio de cada control.
Columnas Evaluación de regla Westgard®	Las reglas Westgard [®] individuales se evalúan como OK si no se ha infringido ninguna regla de control o como x en caso contrario, conforme a la tabla siguiente (ver reglas Westgard a continuación).
Campo Comentarios	Permite escribir comentarios a la evaluación.
Botón Presentación preliminar	Presentación preliminar con el gráfico Levy-Jennings, comprobación de reglas Westgard y comentarios.



Reglas Westgard

Las reglas Westgard® son un modelo QC multiregla:



1:2s	Regla de control para probar si un control de medición ha excedido los límites de control de x + 2SD o x - 2SD.
	Esta regla sirve de alerta para instar a las reglas de control siguientes a una inspección rigurosa de los datos de control.
	Si ninguna medición de control supera la media más 2s o la media menos 2s, entonces el procedimiento de medición se considera " bajo control " aunque se hayan infringido otras reglas.
	Por otro lado, si una medición de control supera la media más 2s o la media menos 2s y se ha infringido alguna otra regla, el procedimiento de medición se considera " fuera de control ".
1:3s	Regla de control para probar si un control de medición ha excedido los límites de control de x + 3SD o x - 3SD .
	Se rechaza un lote si una medición de control supera el límite de control de la media más 3s o la media menos 3s.
2:2s	Regla de control que prueba si dos mediciones de control seguidas exceden el mismo límite de control de x + 2SD o x - 2SD .
	Se rechaza una prueba si 2 mediciones de control seguidas superan el límite de control de la media más 2s o la media menos 2s.
R:4s	Regla de control para probar si el rango o la diferencia entre dos mediciones de control exceden 4SD.
	Se rechaza un lote si una medición de control excede la media más 2s y otra excede la media menos 2s. Los dos resultados de control no necesitan ser consecutivos.
4:1s	Regla de control para probar si cuatro mediciones de control consecutivas exceden el mismo límite de control de x + 1SD o de x - 1SD.
	Se rechaza un lote si 4 mediciones de control consecutivas superan el límite de control de la media más 1s o la media menos 1s.
10:x	Un lote se rechaza si 10 mediciones de control consecutivas caen en el mismo lado de la media. Sin embargo, si uno de estos resultados cae directamente sobre la media, el lote no se rechaza.

7.4.3 Menú de Barra de herramientas: Edición

El menú Edición contiene las siguientes selecciones:

Copiar a Excel

El comando **Copiar a Excel** permite al usuario exportar datos directamente a una hoja de cálculo de Excel de Microsoft.

Copiar en formato ASCII

Seleccione **Copiar en formato ASCII** para copiar al portapapeles los datos contenidos en los pocillos seleccionados del diseño de placa.

Esta función permite transferir los datos de los pocillos a otras aplicaciones.

Pegar desde formato ASCII

En el menú Editar, haga clic en Pegar o pulse Ctrl-V.

El contenido de los datos seleccionados se pegará desde el portapapeles con formato de datos ASCII. Esta función permite transferir datos desde otras aplicaciones de software (por ejemplo, Excel) a los pocillos.

Primero, se deben seleccionar con el ratón en el otro programa los datos relevantes. Después se deben copiar los datos seleccionados para transferirlos al portapapeles. Las filas tienen que estar separadas por avances de línea y las columnas por tabulaciones. Esto se hace automáticamente en Excel cuando se seleccionan varias celdas.

La función **Pegar** pega los datos copiados en formato ASCII del portapapeles a los pocillos. Los datos siempre se insertan empezando por la posición A1. Si los datos no son numéricos, el valor se establece como 0.

En caso de una cinética incompleta con un intervalo definido, se muestra una consulta para preguntar si los datos seleccionados actualmente deberían sobrescribir los datos originales o anexarse a esos datos. El intervalo de tiempo se toma de los parámetros de medición.

Insertar lista de IDs de muestras

Para obtener información detallada, consulte 6.5 Iniciar medición - Insertar lista de IDs de muestras.

Volver a calcular con otro método

Se efectuará un recálculo basándose en los ajustes de un nuevo método seleccionado.

Si se produce un error durante el recálculo, se detendrá el procedimiento.

Haga clic en esta opción para abrir el cuadro de diálogo **Abrir archivo**. En caso de que se haya escrito, en un campo **Notas** adicional se mostrará una descripción de texto del archivo (solamente disponible en **magellan** Tracker). El método se debe seleccionar con doble clic el archivo de método o se debe seleccionar el archivo de método y hacer clic en el botón **Abrir**.

Después de seleccionar el método, se mostrará el cuadro de diálogo **Calculando...** Este cuadro de diálogo solamente es para visualizarlo y no contiene ningún elemento para editar. Se cierra después de que finalice el cálculo.



7.4.4 Menú de Barra de herramientas: Instrumento

Para obtener información detallada acerca de **Movimientos** y **Control de temperatura**, consulte el capítulo 3.1 Opciones del instrumento.

Iniciar medición

Con esta opción, la ejecución de la medición se puede iniciar de nuevo con el método cargado actualmente y los datos actuales se sobrescribirán se ha seleccionado **Sí** en el cuadro de diálogo de **magellan**.

En caso de una ejecución cinética incompleta, se pueden añadir ciclos (no disponible para instrumentos de la serie Infinite).

Para obtener información detallada, consulte 6.5 Iniciar medición con un método predefinido o favorito.

7.4.5 Ventana de diseño de placas

Los datos, que se deben mostrar dentro de un pocillo al abrir un archivo de zona de trabajo, se pueden establecer al definir el método (**Procesamiento** automatizado de datos \rightarrow Ver resultados después de la medición \rightarrow Más...).

En cada pocillo individual, hay tres líneas visibles; de modo predeterminado se mostrarán los siguientes datos:

1ª línea: diseño

2ª línea: información de réplica

3^a línea: en la tercera línea, según los datos seleccionados, se utilizan las normas descritas en el capítulo siguiente (consulte 7.4.6 Caracteres especiales).

7.4.6 Caracteres especiales

"Nada"	No se mostrará ningún dato.
"~" tilde	Se anexa una tilde a los valores manipulado de un pocillo (simulados, editados); por ejemplo: 0,354~
"()" paréntesis	Se aplica paréntesis al valor enmascarado de un pocillo; por ejemplo: (0,354)
"!" signo de admiración	Se anexa un signo de admiración al valor eliminado de un pocillo; por ejemplo: !0,354
"#" almohadilla	Los valores de concentración se marca con una '#' cuando la concentración calculada está fuera del rango de la curva estándar, que significa que el valor de concentración ha sido extrapolado. Por ejemplo: #13,75
"*"asterisco	Un asterisco marca los valores medidos con la opción "Usar regulación de ganancia", que corrige (=reduce) la ganancia.
Se muestra el mensaje de error correspondiente	El valor promedio de un pocillo es inservible (por ejemplo: lámpara tenue, división por cero, logaritmo negativo).
Sin color, la tercera línea está vacía	No hay ninguna fórmula aplicada al pocillo o bien Réplica de pocillo si "promedio s" o "cálculo v" o bien No hay resultado de corte.

Datos de la primera réplica enmascarados o eliminados	Se calcula el valor promedio sin usar datos enmascarados, pero se muestra el valor medio en el pocillo de la primera réplica
"< MÍN"	La concentración calculada está por debajo del estándar inferior.
"MItPt"	Puntos múltiples; no se puede calcular una concentración única.
"> MÁX"	La concentración calculada está por encima del estándar superior.
"NoCalc"	No se encuentran los datos de entrada o error de cálculo; el cálculo no es posible.

Accesos directos

Cuando esté visualizando datos transformados, pulse CTRL-MAYÚS para ver la fórmula definida en la segunda línea mientras pulsa las teclas.

7.4.7 Barra de control: Datos del instrumento

El usuario puede seleccionar datos sin procesar, datos de espectro, etc. y sus estadísticas según la medición realizada.

7.4.8 Barra de control: Datos reducidos

El usuario puede seleccionar datos reducidos, p. ej., datos de diferencia, según la medición realizada.

7.4.9 Barra de control: Datos transformados

El usuario puede seleccionar la transformación requerida para mostrar los resultados calculados y sus estadísticas, presentación coloreada, etc.

7.4.10 Barra de control: Parámetros cinéticos

El usuario puede ver los parámetros de evaluación cinética, como Pendiente Media/Máx., Tiempo hasta iniciación, etc., según la definición de la medición.



Aviso En el cálculo de datos cinéticos, no se tienen en cuenta los datos inservibles (p. ej., los valores de desbordamiento).

Definición de reducción de datos cinéticos específicamente para pocillos

Seleccione **Editar la configuración cinética...** en el menú contextual derecho del ratón cuando tiene un pocillo seleccionado. Se mostrará el cuadro de diálogo Reducción de datos cinéticos (consulte detalles adicionales en la sección 4.3.6 Cinética: Reducción de datos cinéticos).

Para el pocillo seleccionado se pueden definir parámetros particulares que también se pueden copiar a otros pocillos (seleccione **Copiar configuración cinética** en el menú contextual derecho del ratón, haga clic en los pocillos en los que desea copiar la configuración y seleccione **Pegar configuración cinética**...).

Al cambiar la configuración de Reducción de datos cinéticos en la pestaña Método se sobrescribirán las definiciones específicas de los pocillos.

TECAN



Gráfico: Cuadro de diálogo Cinética de enzima

El cuadro de diálogo **Gráfico de cinética de enzima** permite al usuario ver el gráfico de cinética de enzima.

Menú contextual de Gráfico de cinética de enzima

Al hacer clic con el botón secundario en el gráfico se muestra un menú contextual.

Cursor de precisión	Se muestra un cursor de precisión para ayudar a colocar el cursor en puntos particulares sobre el gráfico.
Copiar al portapapeles	Copia el gráfico al portapapeles como un mapa de bits que se puede transferir a cualquier aplicación para Windows utilizando la función pegar.
Guardar como mapa de bits	Guarda el gráfico como un archivo de mapa de bits.
Imprimir	Imprime el gráfico.
Configurar impresora	Muestra el cuadro de diálogo Configurar impresora , en el cual se puede definir la configuración de la impresora.
Zoom al 100%	Si se cambió el factor de zoom seleccionando regiones rectangulares del gráfico, restablece el factor de zoom de la presentación gráfica al 100%.
Datos de cinética de enzima	Abre el cuadro de diálogo Datos de cinética de enzima . El cuadro de diálogo Datos de cinética de enzima muestra el resultado del análisis de cinética de enzima definido.
Propiedades	Seleccione este comando para cambiar las propiedades de presentación del gráfico. Se pueden personalizar el título del gráfico, los estilos del eje, el estilo de la curva, etc.
Ayuda	Abre el cuadro de diálogo de ayuda de magellan.

7.4.11 Barra de control: Concentraciones

Para ver los resultados calculados conforme a la curva estándar, seleccione concentración simple, concentración simple media o promedio:

- Conc. simple (unidad)
 - Concentración de réplica individual.
- Conc. media (unidad)
 Para calcular la concentración se utiliza el valor medio de las réplicas. Si se definen réplicas con distintas diluciones, no estará disponible la concentración media.
- Conc. simple promedio (unidad)
 Se calcula la concentración para cada réplica. A continuación, se calcula un promedio de las concentraciones.
- Hay disponibles concentraciones adicionales (concentración individual, media y promedio).
- Gráfico: Curva estándar Haga clic en esta opción para abrir el cuadro de diálogo Gráfico: Curva estándar
- Intercepciones
- · Valores para transformaciones de concentración, si se han definido
- IC 50, r-IC 50, Gráfico: series de dilución



Gráfico: Cuadro de diálogo Curva estándar

Abra este cuadro de diálogo a través de la barra de control o haciendo clic en **Curva** en la barra de herramientas para mostrar y editar la **Curva estándar**.

Este cuadro de diálogo tiene su propio menú con opciones para guardar la curva estándar, cambiar el tipo de análisis o comparar varias curvas. Ese gráfico se puede exportar como mapa de bits para después poder incorporarlo a documentos de otras aplicaciones.

El cuadro de diálogo **Gráfico: Curva estándar** contiene los siguientes elementos:

Menú Archivo

El menú Archivo contiene los siguientes comandos:

- Nueva se selecciona para crear una nueva curva estándar. Los valores X e Y se pueden definir en un cuadro de diálogo.
 Introduzca valores medidos (valor Y) y de concentración (valor X). Si se selecciona la opción Excluir, el punto correspondiente no se incorporará en el cálculo del gráfico.
 Cuando termine, haga clic en el botón Aceptar.
- Abrir se selecciona para abrir una curva estándar guardada previamente. Se puede abrir una curva estándar guardada en el formato de archivo .std. Al ejecutar este comando, se mostrará el gráfico junto con las curvas estándar ya mostradas.
- Cerrar

Si hay varias curvas abiertas, esta opción permite cerrar curvas individuales. La curva correspondiente se selecciona en la lista y se cierra al hacer clic en el botón **Aceptar**.

• Guardar/Guardar como se selecciona para guardar una curva estándar para realizar evaluaciones adicionales.

Una curva estándar se puede guardar como archivo **.std**. El archivo se puede incluir en el método cambiando a la pestaña **Datos**. Si el usuario selecciona **Guardar** y la curva se había guardado previamente, se guardará sin mostrar ninguna indicación adicional. De lo contrario, se indicará al usuario que introduzca un nombre para la nueva curva.

- **Exportar** se selecciona para exportar una curva estándar como un archivo de mapa de bits.
- **Imprimir...** se selecciona para imprimir la curva estándar. Este comando imprimirá el gráfico actual.
- Configurar impresora...
 Al seleccionar esta opción se muestra el cuadro de diálogo Configurar impresora.
- Salir se selecciona para salir del cuadro de diálogo de curva estándar. Si se ha realizado cualquier modificación, por ejemplo, un cambio en el proceso de interpolación, se realizará un nuevo cálculo.



Aviso

Las curvas estándar agregadas por Nueva o Abrir son solo para realizar comparaciones. Después de cerrar el cuadro de diálogo Curva estándar, esas curvas se eliminarán.

Uso avanzado

En los métodos se puede usar un archivo de curva estándar guardada, que no tenga estándares en el diseño, para calcular concentraciones de muestra. Véase 4.3.8 Concentraciones: Curva estándar.

Menú Editar

El menú Editar contiene los siguientes comandos:

- Seleccione Copiar para copiar el gráfico al portapapeles como un mapa de bits que se puede transferir a cualquier aplicación Windows utilizando la función copiar/pegar.
- Seleccione Propiedades para modificar el tipo de análisis, ejes, título, etc.
- Seleccione Curva estándar para abrir el cuadro de diálogo de curvas estándar. En ese cuadro de diálogo se pueden incluir o excluir los puntos estándar.

Después de seleccionar la curva correspondiente en la lista disponible y hacer clic en el botón **Aceptar**, se mostrarán todos los puntos base en la ventana que aparece. Para activar o excluir un punto, haga clic en la opción **Excluir**. Solo aquellos puntos que no hayan sido marcados como excluidos se usarán al calcular la curva estándar.

Solo se pueden modificar los puntos base obtenidos de datos sin procesar, que han sido introducidos manualmente o que se han cargado adicionalmente. Los puntos base obtenidos, por ejemplo, de transformaciones, solo se pueden excluir.

 Conc. escala: Se muestra el cuadro de diálogo Intervalo de concentración. Elija entre Visualizar todos y Visualizar intervalo. Si se selecciona Visualizar intervalo y se han definido los límites Mín. y Máx., solo se mostrarán en la ventana de diseño de placa los valores de concentración comprendidos en el intervalo especificado.

Menú Ver

El menú Ver contiene los siguientes comandos:

- Seleccione Registro de modificaciones para ver el registro de modificaciones de una curva estándar externa. Esta opción solamente está disponible para magellan Tracker.
- Seleccione **Estadísticas** para mostrar las estadísticas de los valores obtenidos.

Para evaluar si se ha producido alguna desviación o algún cambio en el lector durante un cierto período de tiempo, se pueden comparar series de curvas estándar por medio de los valores estadísticos. Si se cargan varias curvas, se calcularán para cada una de ellas la media, la

Si se cargan varias curvas, se calcularan para cada una de ellas la media, la desviación estándar y el coeficiente de variación.

- Intercepciones... abre el cuadro de diálogo Intercepciones. El cuadro de diálogo Intercepciones muestra los resultados de los valores de la intercepción definida.
- Media de curva estándar muestra la curva estándar promediada sobre los grupos experimentales ejecutados.
 Esta opción solo se puede activar si la ventana presenta varias curvas. Se generará una curva media y se mostrará utilizando los datos de curvas colectivos.

Menú Ayuda

Seleccione Ayuda para abrir el cuadro de diálogo de ayuda de magellan.



Hacer clic en los puntos

Al hacer clic en los puntos, estos se ocultarán/visualizarán dentro de la curva estándar. Después de ocultar un punto, la línea del gráfico se ajusta automáticamente como corresponda y el punto se representa como un símbolo transparente en vez de un símbolo sólido.

Nota: Esto se puede deshacer pulsando Ctrl+Z.

Información sobre herramientas del cursor X, Y

Cuando el cursor no se mueve durante unos instantes, aparece un texto de información sobre herramientas que muestra las coordenadas X e Y de la posición actual del cursor.

Uso avanzado

Si en el gráfico se muestran más de **seis** curvas estándar, solo una pequeña leyenda al lado derecho muestra la etiqueta del gráfico. Para ver las funciones de ajuste con sus parámetros de ajuste de curva respectivos, seleccione **Datos de curva estándar** en el menú contextual (véase a continuación).

Menú contextual de Gráfico de curvas estándar

Al hacer clic con el botón secundario en el gráfico, se muestra un menú contextual.

Cursor de precisión	Se muestra un cursor de precisión para ayudar a colocar el cursor en puntos particulares sobre el gráfico.
Puntos únicos	Muestra valores de réplica en vez de valores medios.
Copiar al portapapeles	Copia el gráfico al portapapeles como un mapa de bits que se puede transferir a cualquier aplicación para Windows utilizando la función pegar.
Guardar como mapa de bits	Guarda el gráfico como un archivo de mapa de bits.
Imprimir	Imprime el gráfico.
Configurar impresora	Muestra el cuadro de diálogo Configurar impresora.
Deshacer cambios	Cancela las acciones más recientes.
Zoom al 100%	Si se cambió el factor de zoom seleccionando regiones rectangulares del gráfico, restablece el factor de zoom al 100% para la presentación gráfica.
Datos de curva estándar	Muestra los parámetros de ajuste de curva de las curvas estándar mostradas.
Intercepciones	Abre el cuadro de diálogo Intercepciones , que muestra los resultados de los valores de la intercepción definida.
Propiedades	Cambia las propiedades de visualización del gráfico. Se pueden personalizar el título del gráfico, los estilos del eje, el estilo de la curva, etc.
Ayuda…	Abre el cuadro de diálogo de ayuda estándar de magellan .



Gráfico: Cuadro de diálogo Series de dilución

El cuadro de diálogo **Gráfico:Series de dilución** muestra el gráfico de dilución con la intercepción especificada.

Menú contextual de Gráfico de series de dilución

Al hacer clic con el botón secundario en el gráfico se muestra un menú contextual.

Cursor de precisión	Se muestra un cursor de precisión para ayudar a colocar el cursor en puntos especiales sobre el gráfico.
Copiar al portapapeles	Copia el gráfico completo al portapapeles como un mapa de bits que se puede transferir a cualquier aplicación Windows utilizando la función pegar.
Guardar como mapa de bits	Guarda el gráfico como archivo de mapa de bits.
Imprimir	Abre el cuadro de diálogo Imprimir , desde el cual se puede imprimir el gráfico.
Configuración de impresora…	Abre el cuadro de diálogo Configurar impresora , en el cual se puede definir la configuración de la impresora.
Zoom al 100%	Si se cambió el factor de zoom seleccionando regiones rectangulares del gráfico, restablece el factor de zoom al 100% para la presentación gráfica.
Datos de series de dilución	Muestra los valores de intercepción y los coeficientes de correlación.
Propiedades	Seleccione este comando para cambiar las propiedades de presentación del gráfico. Se pueden personalizar el título del gráfico, los estilos del eje, el estilo de la curva, etc.
Ayuda	Abre el cuadro de diálogo de ayuda estándar de magellan.

7.4.12 Barra de control: Resultados cualitativos

La definición de corte muestra una perspectiva general de los rangos de corte utilizados actualmente y los límites de corte utilizados.

En la ventana de información de datos de la parte inferior de la pantalla, se muestra la lista de definiciones de corte (grupos experimentales, datos de entrada, rangos de corte, etc.).

El usuario puede ver los resultados de corte.

- Definición de corte Rangos y límites de rangos
- Resultados de corte Resultado de corte para cada pocillo
- Estadística de corte Estadísticas del número de aciertos para cada rango

7.4.13 Barra de control: IDs de muestras

El usuario puede ver los datos de listas de IDs de muestras.



El usuario puede ver el diseño de método, es decir, las posiciones de pocillos, diseño, información de réplicas, parámetros de medición, etc.

TECAN.

7.4.15 Barra de control: Validación QC

La lista de validaciones contiene los grupos experimentales, los datos de entrada para cada grupo de validación, las fórmulas de validación y sus resultados (VERDADERO o FALSO).

Si el resultado es VERDADERO, indica que se cumplieron los criterios de validación para esta placa específica, mientras que FALSO representa un resultado no válido para la placa completa.

7.4.16 Barra de control: Otras opciones

El protocolo de error registra todos los errores que se producen durante la ejecución de un método. Los errores pueden ocurrir durante una medición o durante el cálculo. Revise el protocolo de error antes de utilizar los datos y resultados.

7.4.17 Cuadro de diálogo Escala de colores

Si los valores contenidos dentro de la placa de análisis se presentan en distintos colores, se muestra el cuadro de diálogo **Escala de colores**. Los colores ofrecen una perspectiva general rápida de los resultados de la medición de los pocillos individuales. Los colores, cuando se utilizan, dependen de los ajustes definidos en el juego de colores.

7.4.18 Menú contextual de un pocillo

Si se hace clic con el botón secundario en un pocillo en el diseño de placa (cuando se ha seleccionado la pestaña **Editar método**), se muestra un menú contextual.

Cuadro de diálogo Resumen

El cuadro de diálogo **Resumen** muestra información general de todos los parámetros definidos de un pocillo seleccionado.

Árbol de información	El árbol de información ofrece información general de todos los parámetros basados en pocillos definidos para el pocillo seleccionado. El árbol se rellena según la información disponible, p. ej., si está disponible la información de lista de IDs de muestras, se muestra el ID de muestras, el estado de pipeteado, etc. Si hay disponibles datos sin procesar, se puede mostrar información adicional en función de la configuración de las mediciones y de los instrumentos conectados. Por ejemplo, para las mediciones cinéticas se muestran los puntos de tiempo.
Botones izquierda, derecha, arriba y abajo	Los botones de dirección se pueden usar para seleccionar otro pocillo en la placa mientras se deja el cuadro de diálogo abierto.
Expandir todos / Reducir todos	Haga clic en este botón para mostrar el árbol de información expandido hasta el nivel superior o contraído hasta el primer nivel.
Aceptar	Cierra el cuadro de diálogo Resumen.



Cuadro de diálogo Detalles

El cuadro de diálogo **Detalles** muestra los resultados de múltiples lecturas por pocillo para un pocillo.

Menú Archivo	Guardar y salir se selecciona para guardar todos los cambios y cerrar el cuadro de diálogo.
Menú Editar	 El menú Editar contiene los siguientes comandos: Haga clic en Copiar para copiar valores simples al portapapeles, que entonces se podrán transferir a cualquier aplicación para Windows utilizando la función pegar. Seleccione Volver a calcular para forzar el recálculo de los datos estadísticos del fondo del cuadro de diálogo. Esta opción aparece deshabilitada si se activa Recálculo automático. Si se selecciona Recálculo automático, se calcularán los datos estadísticos después de cada
Menú Ver	 accion de ocultar/visualizar. El menú Ver contiene los siguientes comandos: Se puede seleccionar Valores para mostrar las lecturas como valores. Use esta opción para ver múltiples lecturas por pocillo de forma rápida y aproximada. Se puede seleccionar Gráfico para mostrar los valores leídos como colores. Brillo usa un color con distintos brillos para mostrar los valores utiliza distintos colores para mostrar los valores. Para cambiar la intensidad se muestra una barra de color con un control deslizante.
Menú Ayuda	Haga clic en Ayuda para abrir el cuadro de diálogo de ayuda de magellan .
Hacer clic en los puntos	Al hacer clic en lecturas , estas se ocultarán/visualizarán. Después de ocultar una lectura, su valor se muestra entre paréntesis. Si está habilitada la presentación en color, la lectura se muestra en blanco con un borde circular negro.

Editar diálogo

Utilice este comando para editar los datos sin procesar medidos de un pocillo seleccionado.

Campo de texto Editar datos sin procesar	Se puede introducir un nuevo valor para el pocillo seleccionado. Los datos entonces se muestran con el símbolo ~ (consulte la sección 7.4.6 Caracteres especiales).
Botón Restaurar	Restaura el pocillo a su valor original.
Casilla de verificación Máscara	Excluye del cálculo el valor del pocillo. Entonces los datos se muestran entre paréntesis. Seleccione la casilla de verificación Máscara de nuevo para incluir el valor otra vez.



Botones Izquierda , Derecha, Arriba y Abajo	Los botones de dirección se usan para seleccionar otro pocillo en la placa mientras se deja el cuadro de diálogo abierto.
Aceptar	Cierra el cuadro de diálogo Editar.

Gráfico: Cuadro de diálogo Etiquetas múltiples

El cuadro de diálogo **Gráfico:** El cuadro de diálogo **Etiquetas múltiples** muestra los datos de los pocillos seleccionados. Se muestran los datos sin procesar vs. los ID de los pocillos.

Menú contextual de Gráfico de etiquetas múltiples

Al hacer clic con el botón secundario en el gráfico se muestra un menú contextual.

Cursor de precisión	Se muestra un cursor de precisión para ayudar a colocar el cursor en puntos particulares sobre el gráfico.	
Copiar al portapapeles	Copia el gráfico al portapapeles como un mapa de bits que se puede transferir a cualquier aplicación para Windows utilizando la función pegar.	
Guardar como mapa de bits	Guarda el gráfico como un archivo de mapa de bits.	
Imprimir	Abre el cuadro de diálogo Imprimir desde el cual se puede imprimir el gráfico.	
Configurar impresora	Abre el cuadro de diálogo Configurar impresora , en el cual se puede definir la configuración de la impresora.	
Zoom al 100%	Si se cambió el factor de zoom seleccionando regiones rectangulares del gráfico, restablece el factor de zoom al 100% para la pantalla gráfica.	
Propiedades	Seleccione este comando para cambiar las propiedades de presentación del gráfico. Se pueden personalizar el título del gráfico, los estilos del eje, el estilo de la curva, etc.	
Ayuda	Este botón abre el cuadro de diálogo de ayuda de magellan.	

Gráfico: Cuadro de diálogo Cinética

El cuadro de diálogo **Gráfico:Cinética** muestra los gráficos de mediciones cinéticas de uno o más pocillos seleccionados. La leyenda contiene los parámetros cinéticos calculados.

Haga clic en los puntos para ocultarlos/visualizarlos dentro del gráfico cinético. Después de ocultar un punto, la línea del gráfico se ajusta automáticamente como corresponda y el punto se representa como un símbolo transparente en vez de un símbolo lleno coloreado.

Menú contextual de Gráfico de cinética

Al hacer clic con el botón secundario en el gráfico, se muestra un menú contextual.

Cursor de precisión	Se muestra un cursor de precisión para ayudar a colocar el cursor en puntos particulares sobre el gráfico.
Copiar al portapapeles	Copia el gráfico al portapapeles como un mapa de bits que se puede transferir a cualquier aplicación para Windows utilizando la función pegar.

Guardar como mapa de bits	Guarda el gráfico como mapa de bits.	
Imprimir	Abre el cuadro de diálogo Imprimir desde el cual se puede imprimir el gráfico.	
Configuración de impresora	Abre el cuadro de diálogo Configurar impresora , en el cual se puede definir la configuración de la impresora.	
Zoom al 100%	Si se cambió el factor de zoom seleccionando regiones rectangulares del gráfico, restablece el factor de zoom al 100% para la presentación gráfica.	
Mostrar datos reducidos cinéticos	Los datos cinéticos calculados (pendientes, iniciaciones, mínimos, máximos) se pueden visualizar en el gráfico de curvas cinéticas. Las pendientes se muestran como curvas y las iniciaciones, mínimos y máximos como puntos de intercepción. La presentación se puede activar/desactivar a través del elemento de menú contextual Mostrar datos reducidos cinéticos en el cuadro de diálogo de gráfico cinético. El estado actual de la pantalla se guarda y se reutiliza en futuras presentaciones.	



Curva de temperaturas	Esta opción muestra/oculta una curva que visualiza la progresión de la temperatura durante las mediciones cinéticas (si están disponibles).
Parámetros cinéticos…	Abre el cuadro de diálogo Parámetros cinéticos y ofrece la posibilidad de modificar la configuración cinética de los pocillos seleccionados.
Datos cinéticos	Seleccione esta opción de menú para mostrar una tabla de todos los datos cinéticos calculados de los pocillos seleccionados.
Escala eje Y	Se puede seleccionar un rango para el eje Y.



Propiedades…	Seleccione este botón para cambiar las propiedades de presentación del gráfico. Se pueden personalizar el título del gráfico, los estilos del eje, el estilo de la curva, etc.
Ayuda	Abre el cuadro de diálogo de ayuda de magellan.



Uso avanzado

Es posible establecer distintos parámetros cinéticos para distintos pocillos utilizando el menú contextual para un pocillo en la ventana de diseño de placa o utilizando el menú contextual en el cuadro de diálogo de gráfico cinético. Si se editan los parámetros cinéticos en la pestaña Editar método, los parámetros cinéticos configurados serán idénticos para todos los pocillos.

Gráfico: Cuadro de diálogo Espectros

El cuadro de diálogo **Gráfico: Espectros** muestra el espectro de una medición de exploración.

Menú contextual de Gráfico de espectros 2D

Al hacer clic con el botón secundario en el gráfico, se muestra un menú contextual.

Cursor de precisión	Se muestra un cursor de precisión para ayudar a colocar el cursor en puntos particulares sobre el gráfico.	
Copiar al portapapeles	Copia el gráfico al portapapeles como un mapa de bits que se puede transferir a cualquier aplicación para Windows utilizando la función pegar.	
Guardar como mapa de bits	Guarda el gráfico como un mapa de bits.	
Guardar como JCAMP-DX	Seleccione este comando para guardar el gráfico como un archivo JCAMP-DX 4.24.	
Guardar como ASCII	Seleccione este botón para guardar el gráfico como un archivo ASCII separado por tabuladores.	
Imprimir	Abre el cuadro de diálogo Imprimir , desde el cual se puede imprimir el gráfico.	
Configuración de impresora	Abre el cuadro de diálogo Configurar impresora , en el cual se puede definir la configuración de la impresora.	
Zoom al 100%	Si se cambió el factor de zoom seleccionando regiones rectangulares del gráfico, restablece el factor de zoom al 100% para la presentación gráfica.	
Reducción de datos de espectros	Seleccione este comando para cambiar la configuración de los parámetros de Reducción de datos de espectros .	
Datos reducidos de espectros	Seleccione este comando para mostrar una lista de datos de espectros reducidos. ¡Aviso! Este comando solo está habilitado cuando hay disponibles datos de espectros reducidos de tipo valor.	



Datos como en tabla	Abre un cuadro de diálogo que muestra el espectro en forma de tabla. Esto permite copiar los datos a Excel.
Espectro ocultado	El espectro y los datos reducidos se muestran como ocultos. Ello impide realizar cálculos adicionales para este pocillo.
	Esta opción solo está disponible cuando se muestra el espectro sin procesar.
Mostrar espectro prim.	El espectro seleccionado y el espectro sin procesar se muestran a la vez.
	Esta opción solo está disponible cuando se muestra el espectro reducido.
Propiedades	Seleccione este comando para cambiar las propiedades de presentación del gráfico. Se pueden personalizar el título del gráfico, los estilos del eje, el estilo de la curva, etc.
Ayuda	Abre el cuadro de diálogo de ayuda de magellan.



Aviso

Para exportar los resultados de la exploración, acceda al menú contextual haciendo clic con el botón secundario en el gráfico de espectros. Seleccione Datos como en tabla en el cuadro de diálogo y copie los datos a Excel o guárdelos como archivo ASCII.

7.5 Pestaña Editar método

Haga clic en esta pestaña para cambiar al método utilizado actualmente y su configuración. Cada modificación del método hace que se vuelvan a calcular todos los datos cuando se regrese a la pestaña Evaluar resultados. Estos cambios se pueden guardar en la zona de trabajo. Sin embargo, no se aplicarán al propio archivo de método insertado originalmente.

Consulte el capítulo 4 Asistente Crear o editar método para obtener información adicional.

La pestaña Editar método solamente se muestra si el usuario que esté trabajando en ese momento tiene los derechos apropiados (capítulo 9.6 Derechos del usuario).

7.6 Guardar los Resultados evaluados

Haga clic en **Siguiente** en el cuadro de diálogo **Resultados** para acceder al cuadro de diálogo **Guardar en**:

Cuadro de grupo Guardar en	Campo de texto Nombre de archivo : se mostrará un nombre predeterminado para el archivo, que se puede cambiar.	
	Campo de texto Comentarios del archivo escriba sus notas, si fuera necesario.	
Cuadro de grupo Comentario al registro de modificaciones	Campo de texto Comentarios al registro de modificaciones : escriba sus comentarios en el registro de modificaciones si fuera necesario (solamente disponible para magellan Tracker).	
Botón Guardar:	Guarda el archivo de zona de trabajo	
Exportar	Haga clic en este botón para abrir el cuadro de diálogo Exportación	
Botón Registro de modificaciones:	Se utiliza para ver el Registro de modificaciones del archivo (solamente disponible para magellan Tracker). Para obtener información adicional, consulte el capítulo 4.4 Guardar métodos – Registro de modificaciones.	
Botón Firmas:	Se utiliza para ver las firmas adjuntas al archivo (solamente disponible para magellan Tracker). Para obtener información adicional, consulte el capítulo 4.4 Guardar métodos.	
Firmar esta zona de trabajo ahora	Seleccione Firmar esta zona de trabajo ahora para abrir al asistente Anexar firma (solamente disponible para magellan Tracker).	

Haga clic en el botón **FINALIZAR** para cerrar al asistente **Evaluar resultados**. En caso de que no se hayan guardado los cambios, se preguntará al usuario si desea guardarlos.

Haga clic en el botón Guardar para guardar los datos sin cerrar el asistente.



8. Asistente Anexar firma

8.1 Introducción

A cada registro se le puede aplicar varias firmas, que siempre se incluirán en el informe impreso. Los registros firmados solo los pueden modificar los usuarios con derechos adecuados. Para controlar plenamente la utilización de los métodos se pueden poner límites a los usuarios para que ejecuten exclusivamente métodos firmados. Solo se pueden firmar archivos **Método** y **Espacio de trabajo**.



Aviso Este asistente solamente está disponible con magellan Tracker.

Haga clic en Anexar firma para iniciar el asistente Anexar firma.

Después del cuadro de diálogo de bienvenida, se mostrará el cuadro de diálogo **Seleccionar un archivo**: Seleccione un archivo **Método** o **Espacio de trabajo** que desee firmar.

Cuadro combinado Mostrar	En el cuadro combinado Mostrar se puede modificar la lista de archivos mostrados conforme a la selección. Las selecciones posibles son:
	Todos los archivos
	Archivos sin firma
	Archivos firmados

8.2 Firmar un archivo

Haga clic en Siguiente y se mostrará la ventana Firmar:

rmar		Σ
Firma		
Archivo:	ELISA_Qualitative_M1000.mth	
Significado:	Revisión Acrebación	
	A definir por el usuario:	
Comentario:	for test purposes	
Nombre de usuari Contraseña:	x Admin	
Ayuda Cancelar	<<< Atrás	FIRMAR
		© 2009 Tecan



Confirmación	Solo podrán firmar los usuarios con los derechos requeridos.
Aprobación	Para registros que ya tienen una firma de confirmación, solamente pueden solicitarla los usuarios con derechos apropiados. El mismo usuario no puede firmar la confirmación y la aprobación. Entre la confirmación y la aprobación los archivos no se pueden modificar.

Personalizada Escriba en el área de texto una firma personalizada.

Los ajustes predeterminados del usuario solamente permiten a los administradores anexar firmas a confirmaciones y aprobaciones, así como modificar archivos firmados.

Se pueden añadir comentarios en el cuadro de texto Comentario.

En el campo de texto **Nombre de usuario** se debe escribir el nombre de usuario de la persona que tiene iniciada la sesión en ese momento. En el campo de texto **Contraseña** se debe escribir la contraseña del usuario que tiene iniciada la sesión en ese momento.

Haga clic en **Finalizar** para confirmar la información que ha escrito y firmar el registro.



Aviso

Dependiendo de los procedimientos operativos estándar de la empresa que use el software, esta firma podría tener valor legal. Por lo tanto, de suma importancia que los usuarios mantengan sus contraseñas en secreto.

9. Icono Otras opciones

Haga clic en el icono **Otras opciones** en la ventana del asistente y seleccione entre las siguientes acciones/definiciones:



9.1 Control del instrumento

Consulte el capítulo 3 Control y configuración del instrumento.

9.2 **Procesamiento de archivos**

9.2.1 Guardar en archivador

El cuadro de grupo **Guardar en archivador** ofrece una vista de conjunto y control pleno de la copia de seguridad de los archivos. Contiene una lista de todos los archivos guardados por **magellan** (zonas de trabajo, métodos, listas de IDs de muestras y archivos temporales).



Nombre del archivo	Fechar	Notas		Archivar Seleccionar todo
	20.07.2009.16(28)46	Notas		Seleccional code
21072009-001 wisp	21.07.2009.09:51:19			Recargar
FLISA Sunrise Strin.mth	21.07.2009 10:09:32			Filtro
ELISA Sunrise Evaluat	21.07.2009 10:17:13			Fillero
21072009-003.wsp	21.07.2009 10:17:34			Filtro actual:
ELISA_Sunrise_QC1.mth	21.07.2009 10:37:52			* *
21072009-004.wsp	21.07.2009 10:42:22			
21072009-005.wsp	21.07.2009 10:42:43			
121072009-006.wsp	21.07.2009 11:13:32			
121072009-007.wsp	21.07.2009 11:13:49			
121072009-008.wsp	21.07.2009 11:14:03			
121072009-009.wsp	21.07.2009 11:25:15			
ELISA_Sunrise.mth	21.07.2009 11:41:18			
ELISA_Sunrise_QC.mth	21.07.2009 11:49:51			
21072009-010.wsp	21.07.2009 16:22:31		~	
)		>	
le 64 seleccionados				
overtir	Import	ar	- Loafil	-
			Login	
Convertir de Co	nvertir a Impor	tar datos sin procesar	G	uardar Logfiles

La lista **Nombre de archivo** permite seleccionar archivos para hacer una copia de seguridad de los mismos.

Haga clic en **Archivar** para mover todos los archivos seleccionados a las subcarpetas apropiadas del directorio de copia de seguridad especificado.

Si los archivos tienen el atributo de **solo lectura**, se mostrará un cuadro de aviso antes de eliminarlos.

El cuadro de grupo Guardar en archivador contiene los siguientes elementos:

Directorio Archivador	Defina el directorio que utilizará como Archivador. Haga clic en el botón Examinar para cambiar de directorio.
Lista Nombre de archivo	Contiene todos los archivos guardados por magellan, filtrados según las opciones del botón Filtrar .
Botón Archivar	Haga clic en Archivar para mover todos los archivos seleccionados en la lista Nombre de archivo al directorio de copia de seguridad (función cortar y pegar).
Botón Seleccionar todos	Haga clic en Seleccionar todos para seleccionar todos los archivos que muestra la lista Nombre de archivo .
Botón Recargar	Haga clic en el botón Recargar para anular las opciones actuales de Filtro y actualizar la lista de archivos.
Botón Filtro	 Este botón se utiliza para mostrar solamente archivos específicos, o archivos que tienen una cierta cadena en el nombre de archivo: Se puede utilizar ? como carácter de comodín de un solo carácter. Se puede utilizar * como carácter de comodín de ninguno o varios carácteros



Archivado automático

La función de archivado automático copia a un directorio definido por el usuario los archivos de zonas de trabajo, métodos y muestras que el usuario guarda automática o manualmente.

La opción **Archivado automático** solo está disponible en la versión **magellan Tracker** y solo la pueden habilitar los usuarios con derechos para **Modificar opciones generales**.

tándar
ciones y las
Help

 Haga clic en el botón del explorador de archivos ("...") para examinar el directorio de archivado.



Aviso

Puede suceder que un directorio de archivado se vuelva no válido, p. ej. en una unidad de red. En tal caso, magellan no advertirá explícitamente al usuario, pero añadirá una entrada al archivo de registro de modificaciones del sistema.



Aviso

Tenga en cuenta que los archivos se guardan en el directorio de archivado solo a partir del momento de activar la función Autoarchivar. Los archivos guardados antes de activar la función Autoarchivar no se guardan en el directorio de archivado.



9.2.2 Importar datos sin procesar

La función **Importar de archivo ASCII** permite al usuario procesar en **magellan** archivos de datos ASCII que ya contienen datos de medición y transformaciones adicionales. ES preciso efectuar por adelantado algunas definiciones de formatos de datos para garantizar una extracción de datos correcta. Por ejemplo, los datos individuales deben estar separados de la lista de datos por medio de separadores especificados.

Algunas opciones adicionales que es preciso especificar son: Formato de placa, número de mediciones, intervalo de medición, tipo de orden de datos, orientación y unidad física de los valores.

La importación de datos fallará si el archivo ASCII no contiene tantas entradas como se especifica en el formato de placa.

Una vez que haya definido las opciones requeridas, seleccione el archivo que desee en el campo **selección de archivos** y haga clic en el botón **Abrir** para recuperar los datos del archivo.

El cuadro de diálogo **Abrir** es un cuadro de diálogo estándar de Windows para abrir archivos, con los habituales elementos de navegación entre archivos, un campo de texto para escribir un nombre de archivo y una lista desplegable para seleccionar el tipo de archivo (*.asc).

Además, el cuadro de diálogo **Abrir** contiene los siguientes elementos para establecer opciones ASCII:

Botón de opción Formato	Define el formato de datos previsto, Tabla o Matriz			
Cuadro de grupo Separador	 Define qué carácter se utiliza para separar los campos de datos de la lista o la matriz: El tabulador, si está seleccionado Tabulador. Cualquier otro carácter escrito en el campo de texto Otro. 			
Botón de opción Orientación	Define la orientación de los datos, es decir, horizonta o vertical.			
Parámetros de medición				
Información	 Este cuadro muestra los siguientes parámetros: Formato de placa Unidad de datos Número de mediciones Intervalo de medición 			
Botón Parámetros de medición	Permite seleccionar tipo de medición, tipo de placa, número e intervalo de ciclos cinéticos.			


9.2.3 Convertir a

El cuadro de diálogo **Convertir documentos – Convertir a** permite al usuario convertir archivos de **magellan** desde la versión actual de **magellan** a archivos de versiones anteriores de **magellan**.





Aviso Esta opción solamente está disponible cuando el usuario actual tiene derechos locales de administrador para el sistema Windows.

PRECAUCIÓN

DESPUÉS DE LA CONVERSIÓN, LOS MÉTODOS SE DEBEN VALIDAR PORQUE EL CONTENIDO DE LOS ARCHIVOS CONVERTIDOS PUEDE DIFERIR LIGERAMENTE DEL ARCHIVO ORIGINAL (P. EJ., EL CONTENIDO DEL INFORME IMPRESO, ETC.)

El cuadro de diálogo Convertir documentos contiene los siguientes elementos:

Cuadro de grupo Tipo de documento	Seleccione el tipo de documento que desea convertir
Botón Seleccionar todos	Haga clic en este botón para seleccionar todos los archivos que muestra la lista de archivos
Lista Nombre de archivo	Muestra todos los archivos disponibles para convertir. Para seleccionar un archivo para convertirlo, haga clic en el mismo. Para seleccionar varios archivos, mantenga presionada la tecla de mayúsculas.
Campo de edición Directorio de origen	El campo de edición Directorio de origen contiene el directorio donde se guardan los archivos que se van a convertir. Para cambiar ese directorio, haga clic en el botón a la derecha del campo de texto.
Campo de edición Directorio de destino	El campo de edición Directorio de destino contiene el directorio donde se guardarán los archivos convertidos. Para cambiar ese directorio, haga clic en el botón a la derecha del campo de texto.
Botón Convertir documentos	Haga clic en este botón para iniciar la conversión.
Menú desplegable Versión	Seleccione en ese menú desplegable la versión de magellan a la que se convertirán los archivos.
Barra Progreso	Esta barra muestra el progreso de la conversión.



ADVERTENCIA

LOS ARCHIVOS PUEDEN INCLUIR CARACTERÍSTICAS NO COMPATIBLES CON EL FORMATO SELECCIONADO. EN ESE CASO, ESAS CARACTERÍSTICAS NO ESTARÁN DISPONIBLES.



9.2.4 Convertir de

El cuadro de diálogo Convertir documentos - Convertir de permite convertir archivos magellan de versiones anteriores al formato de archivo actual. También se pueden convertir archivos estándar de magellan a archivos de magellan Tracker.





Solamente se mostrarán archivos para convertir si magellan está conectado al instrumento (en modo de medición o de demostración) para el que se generaron esos archivos.

Aviso







Aviso Deberá convertir los archivos de versiones de magellan anteriores a la 4.0 para poder usarlos.

Aviso Deberá convertir los archivos de magellan Standard para poder usarlos con magellan Tracker.

PRECAUCIÓN

DESPUÉS DE LA CONVERSIÓN, LOS MÉTODOS SE DEBEN VALIDAR PORQUE EL CONTENIDO DE LOS ARCHIVOS CONVERTIDOS PUEDE DIFERIR LIGERAMENTE DEL ARCHIVO ORIGINAL (P. EJ., EL **CONTENIDO DEL INFORME IMPRESO, ETC.)**

El cuadro de diálogo Convertir documentos contiene dos pestañas:

La pestaña magellan 2.x/magellan 3.x contiene los siguientes elementos:

Cuadro de grupo Tipo de documento	Seleccione el tipo de documento que va a convertir, p. ej., un método
Cuadro de grupo Opciones específicas de personalizar nuevo método	Varias opciones globales (en versiones de magellan anteriores a la V4.0) ahora son específicas de los métodos. A fin de establecer estas opciones correctamente en los archivos de método y zona de trabajo, utilice uno de los botones de exportación o el botón "Nombre de zona de trabajo".
Botón Seleccionar todos	Haga clic en este botón para seleccionar todos los archivos que muestra la lista de archivos
Lista Nombre de archivo	Muestra todos los archivos disponibles para convertir. Para seleccionar un archivo para convertirlo, haga clic en el mismo. Para seleccionar varios archivos, mantenga presionada la tecla de mayúsculas. ¡Aviso! Solamente se muestran las zonas de trabajo y los métodos creados con el instrumento conectado
	actualmente.
Campo de edición Directorio de origen	El campo de edición Directorio de origen contiene el directorio donde se guardan los archivos que se van a convertir. Para cambiar ese directorio, haga clic en el botón a la derecha del campo de texto.



Campo de edición Directorio de destino	El campo de edición Directorio de destino contiene el directorio donde se guardarán los archivos convertidos. Para cambiar ese directorio, haga clic en el botón a la derecha del campo de texto. magellan Tracker: este directorio no se puede modificar.
Botón Convertir documentos	Haga clic en este botón para iniciar la conversión.
Barra Progreso	Esta barra muestra el progreso de la conversión.



ADVERTENCIA

LAS PRUEBAS NO CONTIENEN PARÁMETROS DE MEDICIÓN. PARA CONVERTIR UNA PRUEBA EN UN MÉTODO, ES PRECISO DEFINIR LOS PARÁMETROS DE MEDICIÓN. ASEGÚRESE DE QUE LOS PARÁMETROS DE MEDICIÓN SE AJUSTAN A LA PRUEBA SELECCIONADA.

La pestaña **magellan** 4.x / 5.x (disponible solamente con **magellan** Tracker) contiene los mismos elementos de ventana que la pestaña descrita previamente, salvo que las siguientes opciones no son necesarias:

• Probar tipo de documento: magellan 4.x / 5.x no dispone de pruebas.

Opciones específicas de personalizar nuevo método: Estas opciones ya están personalizadas en los archivos **magellan** 4.x / 5.x.

9.2.5 Guardar archivos de registro

Para guardar como archivo de zip los archivos de registro, haga clic en el botón **Guardar archivos de registro...** Entonces podrá escribir un nombre para el archivo zip y guardarlo en un directorio definido. Por si apareciera cualquier error de medición o de estado de **magellan** durante la ejecución, este archivo contiene los datos, los estados (p. ej., desbordamiento o lámpara tenue) y los errores de cálculo, y se puede enviar fácilmente a su departamento de soporte local para solicitar ayuda.



9.3 Opciones

Varias selecciones y ajustes válidos para todos los menús y asistentes de **magellan** se definen en el cuadro de diálogo **Opciones**. Si está activada la administración de usuarios de **magellan**, varios ajustes dependen del usuario. El cuadro de diálogo **Opciones** está dividido en pestañas individuales:

- Pestaña Directorios
- Pestaña Copiar
- Pestaña Vista de placa
- Pestaña Otras opciones

Las opciones solicitadas se deben indicar en las pestañas correspondientes. Haga clic en **Aceptar** para guardar los cambios y cerrar el cuadro de diálogo.

Las opciones que elige el usuario (cada usuario puede modificar esas opciones a su gusto) son:

- Directorios predeterminados (solamente en magellan Standard)
- Opciones de copia al portapapeles y a Excel
- Vista de placa
- Omitir página de bienvenida
- · Empezar con la lista de asistentes o la página de favoritos
- Insertar placa después de comenzar

Las siguientes opciones que dependen del usuario no se pueden definir en el cuadro de diálogo de opciones:

- Impresora seleccionada
- Orientación de impresión.
- Tamaño y fuente de copia impresa.
- Fuente y color de impresión.
- Tipos de identificador predeterminados.
- Dirección de salto para entrada de IDs de muestras.
- Ajustes de Importar datos sin procesar.
- Opciones al exportar predeterminadas, Exportar a archivo ASCII, Exportar a Excel y Nombre de zona de trabajo.
- Tamaño de cuadros de diálogo de gráficos

Las siguientes opciones solamente las puede modificar un usuario con los derechos apropiados. Entonces esas opciones serán válidas para todos los usuarios:

- Idioma
- Permitido Modo de demostración
- Suprimir la advertencia 'no son monótonos' para la curva estándar
- Favoritos
- HUIDs aceptados



Aviso En magellan Standard, el directorio de datos lo puede definir el usuario. En magellan Tracker el directorio de datos es un ajuste general, válido para todos los usuarios.



9.3.1 Directorios predeterminados de datos

Esta pestaña define las rutas de acceso predeterminadas a los directorios en los que se guardan los distintos tipos de archivos utilizados en **magellan**. Estas rutas de acceso predefinidas se configuran al instalar el software **magellan** y se pueden cambiar en los campos de datos correspondientes de la pestaña **Directorios**.

C:\Users\Pi				
	ublic\Docume	ents\Tecan\l	Magellan\wsp	
Método:				
C:\Users\Pi	ublic\Docume	ents\Tecan\l	Magellan \mth	
Lista de IDs	de muestra:			
C:\Users\Pi	ublic/Docume	ents\Tecan\l	Magellan\smp	
Exportación:				
C:\Users\Pi	ublic/Docume	ents\Tecan\l	Magellan \asc	
Curva estáno	dar:		a a a	-
C:\Users\Pi	ublic \Docume	ents\lecan\l	Vlagelian \wsp	
Bitmaps:			Maaallaa	
C. \Users \Ft	JDIIC (DOCUME	ints v recari v	viagelian	
	Cambia	r directorio pr	edeterminado de	datos

La pestaña Directorios contiene los siguientes elementos:

Campos de texto Directorios predeterminados	Se muestra un campo de texto con la ruta predeterminada correspondiente a cada uno de los siguientes tipos de archivo:
	Zona de trabajo
	Método
	Lista de IDs de muestra
	• Exportación: para datos de medición exportados a archivos ASCII.
	 Curva estándar: la ruta para las curvas estándar es siempre la misma que la de las zonas de trabajo.
	 Bitmaps: para mapas de bits creados en los cuadros de diálogo de gráficos.
Botones 🛄	Haga clic en los botones que hay junto a los campos de texto para abrir un cuadro de diálogo estándar de Windows donde se puede seleccionar una ruta de acceso distinta para los archivos generados.



Botón Cambiar directorio predeterminado de datos Solo disponible en **magellan** Tracker. Haga clic en este botón para cambiar la ruta predeterminada para guardar los archivos de zonas de trabajo, métodos, listas de IDs de muestras y curvas estándar. Para modificar los directorios de datos predeterminados se necesitan derechos de administrador y de administrador del sistema de **magellan**.



ADVERTENCIA

ES POSIBLE QUE EL SOFTWARE NO PUEDA ASIGNAR LOS DERECHOS CORRECTOS DEL SISTEMA DE ARCHIVOS A UNA RUTA DE DATOS PREDETERMINADA DEFINIDA POR UN USUARIO. ESA OPERACIÓN DEBE REALIZARLA MANUALMENTE EL ADMINISTRADOR DEL SISTEMA. NO SE DEBE PERMITIR BORRAR O MODIFICAR ARCHIVOS EN LOS DIRECTORIOS PREDETERMINADOS.



Aviso En Magellan Tracker solamente se pueden modificar las rutas para Exportación y Bitmaps.

9.3.2 Opciones de Copiar/Exportar

En un documento de zona de trabajo, los comandos **Editar – Copiar** y **Editar – Copiar a Excel** permiten al usuario exportar datos de mediciones al portapapeles o directamente a una hoja de cálculo Excel. En la pestaña **Copiar** se especifica de qué modo se procesa una placa en caso de transferencia de datos.



150



Cuadro de grupo Formato de matriz	El usuario también tiene las siguientes posibilidades de definir cómo desea que se muestren los datos.
	 Botón de opción Filas en orden horizontal Las filas de datos de la microplaca se copiarán en orden horizontal.
	 Botón de opción Filas en orden vertical Las filas de datos de la microplaca se copiarán en orden vertical. Esto dará como resultado una matriz transpuesta.
	 Si está seleccionada la casilla de verificación Añadir etiquetas de fila/columna, se añade la información de fila y columna como en XFluor cuando se copian los datos.
Cuadro de grupo Opciones de copia	Se debe seleccionar un formato para el punto decimal de los números de los datos transferidos:
a portapapeles	 Botón de opción Punto (x.xxx) Ejemplo: 7.893
	 Botón de opción Coma (x,xxx) Ejemplo: 7,893
Cuadro de grupo Opciones de copia	Los botones de opción permiten definir dónde situar los datos transferidos dentro de Excel.
a Excel	 Botón de opción Nuevo libro de trabajo Los datos transferidos se guardarán en la primera hoja de cálculo de un nuevo libro de trabajo. Siempre se creará un nuevo archivo, independientemente de si Excel está abierto o no.
	 Botón de opción Nueva hoja de cálculo Los datos transferidos se ubicarán en una nueva hoja de cálculo dentro de un libro de trabajo Excel abierto v activo.
	Al seleccionar esta opción, si Excel no está abierto se abrirá, se creará un nuevo libro y los datos solicitados se insertarán en la primera hoja de cálculo.
	 Botón de opción Insertar en hoja de cálculo en la celda y campo de texto de coordenadas de celda Los datos transferidos se ubicarán en la celda indicada (la celda predeterminada es A1) de una hoja de cálculo Excel abierta y activa. Al seleccionar esta opción, si Excel no está abierto se abrirá, se creará un nuevo libro y los datos solicitados se insertarán en la primera hoja de cálculo.
	 Botón de opción Anexar a la hoja de cálculo actual Esta opción simplemente anexará los datos a la hoja de cálculo actual. Al seleccionar esta opción, si Excel no está abierto se abrirá, se creará un nuevo libro y los datos solicitados se insertarán en la primera hoja de cálculo.

La pestaña Copiar contiene los siguientes elementos:



9.3.3 Pestaña Vista de placa

En la pestaña **Vista de placa** se puede modificar el aspecto de la vista de placas. Se podrán cambiar los colores de las diferentes áreas mediante la selección del color deseado en la lista desplegable. El modo de selección para marcar pocillos también se puede definir aquí. También se puede ajustar el tamaño de fuente.

lirectorios	Copiar	Vista de plac	a Otras	opciones		
						^
	0	1	2	3	4	1
	A					_
	В					
	C					Γ
<		-				~
Fondo		•	Sombr	ra clara		-
Fondo de p Superficie	olaca [Sombr	ra oscura		 +
placa			1		0	Unión

La pestaña Vista de placa contiene los siguientes elementos:

Botón Fondo	Para elegir el color de fondo de la ventana Visualizar placa .
Botón Fondo de placa	Para elegir el color del marco de la placa.
Botón Superficie de placa	Para elegir el color de la superficie de la placa.
Lista desplegable Tamaño fuente	Para elegir la altura de letras del texto en los pocillos:PequeñoMedioGrande
Botón Sombra clara	Para elegir el sombreado claro del marco de la placa.
Botón Sombra oscura	Para elegir el sombreado oscuro del marco de la placa.



Botones de opciónAquí se establece el modo de selección para marcarModo de selecciónpocillos con el ratón:

- **Unión** permite seleccionar los pocillos solo cuando se ha marcado el área completa de estos con el ratón.
- **Cruce** permite seleccionar los pocillos solo cuando se ha marcado una parte del área de estos con el ratón.

9.3.4 Otras opciones

pciones			2
Directorios	Copiar	Vista de placa	Otras opciones
Asistente ⊠Omitir bienvo	página enida	de	Iniciar al comienzo Cuadro de diálogo Lista de asistentes Cuadro de diálogo Favoritos
Permitie Meter p Suprim Establece	do Modo blaca de ir la adv r siguier	o de demostraci espués de come ertencia 'no son nte valor del cor	ón nzar monótonos' para la curva estándar itador
Idioma Seleccion	nar: E	spañol	reinicio diario:
Entomos i Aceptar a	regulado archivos	os se ordenadores	s (HUID): Añadir HUIDs
Almacena	miento a ar almac	automático enamiento	
Cambiar o	contrase	ña Regi	stro de modificaciones de opciones y onfiguraciones predeterminadas
		ОК	Cancel Help

La pestaña Otras opciones contiene los siguientes elementos:

Cuadro de grupo Asistente	Cuando se selecciona la casilla de verificación Omitir página de bienvenida , al abrir un asistente no se muestra la página de bienvenida del asistente. El asistente comienza en su segunda ventana.
Cuadro de grupo Iniciar al comienzo	Al seleccionar la casilla de verificación Cuadro de diálogo Favoritos se establece que en lugar de la <i>Lista</i> <i>de asistentes</i> se mostrará el <i>Cuadro de diálogo Favoritos</i> después de iniciarse el software magellan .
Cuadro de grupo	Casilla de verificación Permitido Modo de demostración : Seleccione esta opción para permitir la conexión a un instrumento simulado. Esta opción resulta muy práctica si se desea evaluar datos ya medidos. Entonces no se necesitará ningún instrumento.



	 Casilla de verificación Meter placa después de comenzar: Si se selecciona esta casilla de verificación y se conecta un instrumento después de iniciarse, el portaplacas se introducirá automáticamente.
	 Casilla de verificación Suprimir advertencia "no son monótonos": Elija esta opción para suprimir la advertencia "no son monótonos". Esta opción es útil si en una curva estándar se toman varios puntos en un área plana y es muy probable que no ofrezca monotonía (es decir la parte plana de una curva logarítmica u otra curva donde se aproxima un límite). ¡Aviso! Esta opción está deshabilitada en magellan Tracker.
Cuadro de grupo Establecer siguiente valor del contador	 Se ofrecen dos opciones: El control de entrada numérica Contador total permite establecer el contador a cualquier valor. Muestra el número asignado al siguiente archivo de zona de trabajo creado. El control de entrada numérica Contador con reinicio diario permite establecer el contador a cualquier valor. Muestra el número actual de los archivos de zona de trabajo creados durante el día.
Cuadro de grupo Idioma	El campo de texto muestra el idioma aplicado. Se puede elegir otro idioma en la lista desplegable. Véase Añadir a continuación.
Botón Añadir	Seleccione otro idioma y haga clic en el botón Añadir para reiniciar el programa con el idioma elegido.
Botón Cambiar contraseña	Haga clic en este botón para abrir un cuadro de diálogo donde el usuario puede cambiar su contraseña (solo disponible con Administración de usuarios de magellan activada).
Botón Añadir HUIDs	Se puede definir una lista adicional de HUIDs que permite al usuario abrir archivos creados en otro ordenador (solo disponible en magellan Tracker).
Archivado automático	Seleccione esta función y especifique un directorio definido por el usuario para copiar todos los archivos de zona de trabajo, método e ID de muestras que el usuario guarda manual o automáticamente en un directorio de archivado definido por el usuario.
Botón Registro de modificaciones de opciones y configuraciones predeterminadas	 Haga clic en este botón para ver el registro de modificaciones en opciones y configuración predeterminada. El registro de modificaciones se puede guardar en un archivo ASCII (solo disponible en magellan Tracker). Haga clic en Guardar como archivo para guardar esta información como archivo .log. o .txt. Ese texto también se puede copiar en un procesador de textos para imprimirlo.



9.4 Administración de usuarios (magellan Tracker)

En este documento se mencionan tres tipos de administradores:

El **Administrador del sistema** se encarga de los cambios necesarios en el sistema operativo del ordenador.

El **Administrador** es el responsable de los derechos del usuario en el software **magellan**.

El **Administrador** de **magellan** tiene acceso a todos los componentes del software **magellan**, pero no puede cambiar los derechos del usuario.

magellan ofrece protección mediante contraseña para evitar el uso indebido del software y para restringir el acceso a ciertas partes del software según los derechos de usuario.



 Image: Registro de modificaciones
 Image: Opciones
 Image: Resumen
 Image: Opciones
 Image: Opcione

Roles	La lista Roles muestra todos los grupos de usuarios (niveles de usuarios) existentes.
Agregar rol	Permite crear un nuevo grupo con los derechos correspondientes.

contiene los siguientes elementos:

🛵 Cambiar contraseña



Modificar rol	Permite modificar los derechos de un grupo ya existente.
Usuarios	Esta lista muestra los usuarios existentes y los grupos a los que pertenecen.
Añadir usuario	Permite crear un nuevo usuario.
Modificar usuario	Permite cambiar el nombre completo, el nivel de usuario (grupo) o la contraseña de un usuario ya existente.
Bloquear usuario	Permite habilitar o deshabilitar una cuenta de usuario. Este botón solamente será visible si hay un nombre de usuario seleccionado.
Registro de modificaciones	El servidor de gestión de usuarios escribe en el registro de modificaciones de gestión de usuarios todos los cambios realizados en la base de datos de administración de usuarios (p. ej., creación de grupos/usuarios, cambio de derechos, cambio de opciones, etc.).
Opciones	Permite modificar las opciones generales de inicio de sesión, contraseña o correo electrónico.
Resumen	Muestra una descripción detallada de todos los usuarios y grupos de usuarios existentes como texto sin formato

9.4.1 Añadir/modificar usuario (magellan Tracker)

Haga clic en **Añadir usuario** en el cuadro de diálogo **Administrar usuarios y roles**. Se mostrará el cuadro de diálogo **Crear usuario**:

🥜 Crear usuario					
Crear usuario Escriba el nombre de usuario, la contraseña y el nombre completo del usuario que desea crear.					
Nom <u>b</u> re de usuario:	ToEx				
No <u>m</u> bre completo:	Tom Example				
Contraseña:	***** ?				
C <u>o</u> nfirmar contraseña:	****				
	<u>A</u> ceptar Ca <u>n</u> celar				

El cuadro de diálogo Crear usuario contiene los siguientes elementos:

Nombre de usuario	Para crear un nuevo usuario escriba un nuevo nombre de usuario. Este nombre no se podrá modificar más tarde.
Nombre completo	Escriba el nombre completo del usuario. Podrá cambiarlo más tarde.
Contraseña	Escriba la contraseña inicial. Esta contraseña se deberá cambiar en el próximo inicio de sesión.



Haga clic en Aceptar y se mostrará el cuadro de diálogo Modificar usuario:

💄 Modificar usuar	io				
Actualizar usuario Actualiza el usuario: cambia su nombre completo, contraseña y rol asociado. El nombre de usuario no se puede cambiar.					
Cambiar imagen Eliminar imagen	N <u>o</u> mbre de usuario: <u>N</u> ombre <u>R</u> oles:	ToEx Tom Example Administrators Magellan_Operator Magellan_Application_Specialist Magellan_Administrator Users			
		∫ <u>x ⊆</u> ambiar contraseña <u>A</u> ceptar Canc <u>e</u> lar			

El cuadro de diálogo Modificar usuario contiene los siguientes elementos:

Nombre completo	Escriba el nombre completo del usuario.		
Roles	Seleccione uno de los grupos de usuarios existentes. Al usuario se le asignarán los derechos asociados con ese grupo.		
Cambiar imagen	Se puede agregar o cambiar la foto del usuario.		
Eliminar imagen	Borra la foto del usuario.		
Cambiar contraseña	Se puede definir una nueva contraseña de usuario. Tenga en cuenta que el administrador no necesitará introducir la contraseña antigua.		

Para modificar un usuario, haga clic en **Modificar usuario** en el cuadro de diálogo **Administrar usuarios y roles**. Acceda al cuadro de diálogo **Modificar usuario** y prosiga como se ha descrito.

9.4.2 Agregar/modificar grupo

Haga clic en **Agregar rol** en el cuadro de diálogo **Administrar usuarios y roles**. Se mostrará el cuadro de diálogo **Crear nuevo rol**: Defina un **Nombre de rol** único. Tenga en cuenta que ese nombre no se podrá modificar después. Haga clic en **Aceptar** para cerrar el cuadro de diálogo **Agregar rol**. Abra el cuadro de diálogo **Modificar rol**:

🐴 Modificar rol		×		
Actualizar rol Para actualizar el rol, modifique las asociaciones con los derechos de aplicación. El nombre del rol no se puede modificar.				
<u>N</u> ombre de rol:	Users			
Aplicación:	Magellan	*		
<u>D</u> erechos:	 Create sample id lists Edit sample id lists Sign sample id lists Modify signed sample id lists Create methods Edit methods Sign methods Modify signed methods 	•		
	<u>A</u> ceptar Canc <u>e</u> lar	·		

El cuadro de diálogo Modificar rol contiene los siguientes elementos:

Nombre del rol	El nombre del rol se define en el cuadro de diálogo Agregar rol.
Aplicación	Componente del software de Tecan, p. ej. Magellan.
Derechos	El rol representa el nivel de un usuario determinado. Se pueden seleccionar los derechos de usuario correspondientes.



9.4.3 Registro de modificaciones de administración de usuarios

Haga clic en **Registro de modificaciones** en el cuadro de diálogo **Administrar usuarios y roles**. Se mostrará el cuadro de diálogo **Registro de modificaciones**:

🗉 🗀 C:\Prog	ramData\Tecan	NTecan User M	anagement\v1.	I \Audit \	
	uments and Sett 118-02-10 09-28-28	tings\All Users\ -380.log	Application Data	a\Tecan\Tecan l	Jser Managem
- 20 20	18-02-10 09-28-28	-661.log			
	18-02-10 09-28-29 18-02-10 09-28-58	-708.log -292.log			
1					
6	🔰 = checksum c	of file is correct	🛞 = checksun	n of file is incorrect	
\$ \$	🕅 = checksum c	of file is correct	장 = checksun	n of file is incorrect	_
this selected file	> = checksum o was modified	of file is correct	장 = checksun	n of file is incorrect	
This selected file	= checksum o	of file is correct	🛞 = checksun	n of file is incorrect	
This selected file by on	> = checksum o	of file is correct	😿 = checksun	n of file is incorrect	
This selected file by on	> = checksum o	of file is correct	😿 = checksun	n of file is incorrect	

El servidor de gestión de usuarios escribe en el registro de modificaciones de gestión de usuarios todos los cambios realizados en la base de datos de administración de usuarios.

Para ver un archivo de registro, selecciónelo en la lista (seleccione la casilla de verificación) y haga clic en el botón **Imprimir**. El archivo se abrirá en el visor de PDF y entonces se podrá imprimir. La marca de verificación verde indica que los archivos tienen una suma de control correcta y el aspa roja indica que dicha suma es incorrecta.



Aviso

Tenga en cuenta que el archivo de registro actual siempre tiene una aspa roja porque el archivo aún se estará modificando mientras el software esté abierto.

Derechos del usuario

En los derechos de usuario se selecciona el nivel de seguridad entre los tres existentes. El nivel más alto es **Administrador de Magellan**, que tiene acceso a todas las funciones del programa. Los niveles **Especialista en aplicaciones** y **Operador** tienen más limitaciones.

magellan Standard: se asignan los derechos predeterminados como se detalla a continuación.

magellan Tracker: un Administrador puede asignar o retirar cada uno de los derechos y de modo predeterminado se asignan como se detalla a continuación. El conjunto predeterminado de derechos está basado en las condiciones típicas de laboratorio.





PRECAUCIÓN

LOS ADMINISTRADORES TIENEN LA RESPONSABILIDAD DE VERIFICAR QUE SE ACATAN LAS LEYES APLICABLES Y QUE LOS AJUSTES (Y TODAS LAS MODIFICACIONES) SE CORRESPONDEN CON LAS CONDICIONES DE LABORATORIO DEFINIDAS EN LOS PROCEDIMIENTOS OPERATIVOS ESTÁNDAR (SOP) DE LOS USUARIOS.



Aviso En el caso de aplicaciones de diagnosis clínica, la autoridad competente debe validar todos los métodos para garantizar la seguridad de los resultados. Un método se considerará validado cuando está firmado por la autoridad competente. Dicha autoridad es totalmente responsable de los resultados obtenidos.

9.4.4 Opciones de Administración de usuarios

Haga clic en **Opciones** en el cuadro de diálogo **Administrar usuarios y roles.** Se mostrará el cuadro de diálogo **Opciones de edición**.

Opciones de inicio de sesión

Opciones de ed	ción			×		
Iniciar sesión	Contraseña	Correo electrónico				
Bloquearla	aplicación de	espués de estos minutos de inactivida	d:			
15		-				
Salir de la de este nú	aplicación y bl mero de intent	loque <mark>a</mark> r la cuenta después tos fallidos de inicio de sesión:				
3		÷				
Bloquear aplicación nicios de	La (ur usu De	aplicación se bloquea de n minuto a 36500 h). Par uario debe escribir de nu spués de varios intentos	OK Cancelar espués de un tier a desbloquear la uevo su contraseñ s de inicio de sesi	npo aplic ía. ón fa	especifi cación, d	cado ∍l
Bloquear aplicación nicios de sesión allidos	La (ur usu De 100 env adu	aplicación se bloquea de n minuto a 36500 h). Par uario debe escribir de nu spués de varios intentos 0), la cuenta del usuario vía un mensaje de correc ministrador del sistema.	OK Cancelar espués de un tier a desbloquear la uevo su contraseñ s de inicio de sesi se bloquea y opo o electrónico de a	npo aplic ía. ón fa ciona alerta	especifi cación, o allidos (almente a al	cado el 1 – se





PRECAUCIÓN

SI SE OLVIDA LA CONTRASEÑA DEL ADMINISTRADOR DE USUARIOS, NO SE PODRÁN CAMBIAR LOS AJUSTES DE LA ADMINISTRACIÓN DE USUARIOS. EN ESE CASO, PRIMERO DESINSTALE MAGELLAN Y LUEGO EFECTÚE UNA NUEVA INSTALACIÓN (CONSULTE LA SECCIÓN 1.5 PROCEDIMIENTO DE INSTALACIÓN DEL SOFTWARE).



Aviso Al desinstalar magellan, no se perderá NINGÚN archivo de magellan existente. Todos los archivos previos se podrán encontrar en una carpeta de respaldo generada durante el nuevo procedimiento de instalación.

Opciones de contraseña

niciar sesión	Contraseña	Correo electrónico		
Longitud m	ínima requerio	la de la contraseña:	Permitir contras	eña vacía inistrador)
Número m	ínimo requerido	o de caracteres no alfani	iméricos:	
0	100	-		
Número m	ínimo requerido	o de dígitos:		
0		+		
Número m	ínimo de camb	ios <mark>de contraseña para p</mark>	oder reutilizar una:	
3		+		
La contras	eña caduca d	espués de este número c	le días:	
90				
			OK	Cancelar

Las contraseñas deben contener un número mínimo especificado de caracteres. Se puede definir que las contraseñas contengan caracteres no alfanuméricos y/o dígitos. No se pueden volver a usar las contraseñas ya usadas.

Longitud mínima de la contraseña:

Especifique el número requerido de caracteres.

Número mínimo requerido de caracteres no alfanuméricos:

Especifique el número requerido de caracteres no alfanuméricos.

Número mínimo necesario de dígitos:

Especifique el número requerido de dígitos.

Número mínimo de cambios de contraseña para poder reutilizar una:

Especifique el número de cambios de contraseña.

La contraseña caducará después de este número de días:

Las contraseñas de usuario caducan después de un número configurable de días. Entonces el usuario deberá seleccionar otra contraseña. Especifique el número requerido de días (entre uno y 36500 días).



Permitir contraseña vacía:

Seleccione esta casilla de verificación si desea que se permita que un usuario tenga la contraseña en blanco (excepto para el administrador).

Opciones de correo electrónico

Notificar por correo electrónico	
Enviar por correo electrónico a:	
Enviar por SMTP Servidor	
Puerto	:
Protocolo de seguridad:	○ SSL ○ TLS
Remite	:
Autenticación	: []
Usuario	6

Se notificará al administrador si se producen posibles ataques a la seguridad del sistema (se ha bloqueado una cuenta de usuario debido a varios inicios de sesión incorrectos).

Se puede enviar un mensaje de correo electrónico a la dirección de correo electrónico escrita:

• Servidor SMTP: se debe escribir la dirección IP del servidor SMTP (p. ej., SEUATEXCH01.eu.tecan.net)

Haga clic en **Correo de prueba** para enviar un mensaje de prueba para verificar que los ajustes son correctos.



Aviso La notificación a través de correo electrónico requiere un servidor SMTP sin autenticación.

9.4.5 Resumen de Administración de usuarios

Para obtener una descripción detallada en forma de texto de todos los usuarios y grupos de usuarios existentes, en el cuadro de diálogo **Administrar usuarios y roles,** haga clic en **Resumen**.

Ese texto se podrá copiar a un programa procesador de textos para imprimirlo.



9.5 Administración de usuarios (magellan Standard)

El software **magellan** dispone de una protección por contraseña para evitar un mal uso del software y el acceso por parte de usuarios no autorizados.

A	dministración de usuarios		
	Nombre de usuario Admin (Admin) APSP (Application Specialist)	Derechos Administrador_Ma Operador_Magellan	Añadir usuario Modificar Desactivar
	Protección por contraseña (activ	ar/desactivar)	Iniciar sesión
		Aceptar Cancelar	Ayuda

Los perfiles de usuario se pueden definir con tres niveles de seguridad distintos. El nivel de seguridad más alto dentro del software es el de **Administrador** de **magellan**. Este nivel otorga acceso a todas las funciones del programa y de edición de datos. Para los niveles de **Especialista en aplicaciones** y **Operador**, las opciones son cada vez más restringidas.

Solamente el **Administrador de usuarios** tiene la capacidad de habilitar o deshabilitar la protección por contraseña del software, definir nuevos usuarios o eliminar los antiguos. Si los usuarios han iniciado sesión con su nombre de usuario, el sistema solamente dará acceso a las opciones conforme a su definición de derechos de usuario.

El cuadro de diálogo **Administración de usuarios** contiene los siguientes elementos:

Botón Inicio de sesión	Se mostrará el cuadro de diálogo Inicio de sesión . Los usuarios pueden iniciar sesión con su nombre de usuario.
Lista de usuarios	En esta lista se muestran todos los usuarios definidos en ese momento con su Nombre de usuario y Derechos . Aquí se puede seleccionar cualquier usuario para modificarlo o eliminarlo.
Botón Añadir usuario	El botón Añadir usuario abre el cuadro de diálogo Derechos del usuario , en el que se deberá escribir el ID de usuario, el nombre y la contraseña para definir un nuevo usuario.
Botón Modificar	El botón Modificar abre el cuadro de diálogo Derechos del usuario para poder modificar el usuario seleccionado en la lista de usuarios.
Botón Desactivar	El botón Desactivar deshabilitará el usuario seleccionado en la lista de usuarios actual.
Casilla de verificación Protección por contraseña (activar/desactivar)	Esta casilla de verificación permite habilitar o deshabilitar la función de protección por contraseña del software. Si está habilitada, se mostrará el cuadro de diálogo Inicio de sesión al iniciarse magellan.

9.5.1 Añadir/modificar usuario (magellan Standard)

Este cuadro de diálogo permite añadir o modificar usuarios y asignarlos a un Grupo de usuarios. Para conocer los derechos de los distintos grupos de usuarios, consulte el capítulo 9.6 Derechos del usuario.

El cuadro de diálogo Derechos del usuario contiene los siguientes elementos:

Campo de texto Nombre de usuario	Escriba el nombre completo del usuario, tal como desea mostrarlo
Campo de texto ID de usuario	Defina el ID de usuario que el usuario tendrá que escribir para iniciar sesión
Botones de opción Derechos del usuario	 Defina el nivel de acceso a las funciones del software disponibles para el usuario: Operador Especialista en aplicaciones Administrador
Campo de texto Contraseña	Defina la contraseña del usuario (al menos cinco caracteres)
Campo de texto Confirmar contraseña	Escriba de nuevo la contraseña. Por seguridad, en este campo los caracteres de la contraseña se muestran como un símbolo.

Haga clic en el botón Aceptar para guardar y cerrar el cuadro de diálogo.

Haga clic en el botón **Cancelar** para cerrar el cuadro de diálogo sin guardar los ajustes que ha escrito.



Si no se ha definido todavía ningún usuario y se va a usar Administración de usuarios, primero se debe definir un administrador. Por lo tanto, al hacer clic en el botón Inicio de sesión, se mostrará un mensaje corto. Después se mostrará el cuadro de diálogo Derechos del usuario para poder escribir la información de usuario del Administrador (consulte el capítulo 2.3 Iniciar magellan/ Inicio de sesión).

9.5.2 Inicio de sesión

Este botón solamente estará habilitado si está activada la Administración de usuarios, pero no se ha seleccionado **Protección por contraseña**. Por lo tanto, en el inicio no se muestra el cuadro de diálogo **Inicio de sesión**. A fin de activar la administración de usuarios protegida por contraseña, haga clic en el botón **Inicio de sesión**.

9.5.3 Cambiar usuario

Para cambiar el usuario activo, haga clic en el icono **Cambiar usuario** en la ventana de asistentes. Se mostrará el cuadro de diálogo **Inicio de sesión**.



9.6 Derechos del usuario

Existen tres niveles de seguridad de derechos de usuario. El nivel más alto es **Administrador**, que tiene acceso a todas las funciones del programa. Los siguientes niveles, **Especialista en aplicaciones** y **Operador** tienen mayores limitaciones.

magellan Standard: se asignan los derechos predeterminados como se detalla a continuación.

magellan Tracker: un administrador de usuarios puede asignar o retirar cada derecho y de modo predeterminado se asignan como se detalla a continuación. El conjunto predeterminado de derechos está basado en las condiciones típicas de un laboratorio.





LOS ADMINISTRADORES DE USUARIOS TIENEN LA RESPONSABILIDAD DE GARANTIZAR QUE LOS AJUSTES (Y LAS MODIFICACIONES) SE CORRESPONDAN CON LAS CONDICIONES DE LABORATORIO DEFINIDAS EN LOS PROCEDIMIENTOS OPERATIVOS ESTÁNDAR (SOP) DE USUARIOS Y QUE CUMPLAN LAS LEYES APLICABLES.



Aviso

En el caso de aplicaciones de diagnosis clínica, la autoridad competente debe validar todos los métodos para garantizar la seguridad de los resultados. Un método se considerará validado si está firmado por la autoridad competente. Dicha autoridad es totalmente responsable de los resultados obtenidos.

Se han resumido los derechos del usuario en la siguiente tabla, y son válidos para ambas versiones de **magellan**, es decir. **magellan Tracker** y **magellan Standard**, a menos que se indique lo contrario:

9. Icono Otras opciones



Derechos del usuario	Administrador de magellan	Especialista en aplicaciones	Operador de magellan
Lista de IDs de muestras		ue magenan	
Crear listas de IDs de muestras			
Crear una nueva lista de IDs de muestras o guardar una lista ya existente con un nuevo nombre.	Sí	Sí	Sí
Editar listas de IDs de muestras			
Editar una lista de IDs de muestras ya existente y guardarla con el mismo nombre.	Sí	Sí	No
Firmar listas de IDs de muestras*			
Firmar la revisión/aprobación de una lista de IDs de muestras ya existente.	Res	ervado para uso fut	uro.
Modificar listas de IDs de muestras firmadas*			
Editar o cambiar IDs de muestras en una lista de IDs de muestras firmada.	Res	ervado para uso fut	uro.
Métodos			
Crear métodos			
Crear un nuevo método o guardar con un nuevo nombre un método existente.	Sí	Sí	No
Editar métodos			
Editar un método existente y guardarlo con el mismo nombre.	Sí	Sí	No
Firmar métodos*			
Firmar un método existente para revisión/aprobación.			
Nota: El derecho Firmar métodos por sí solo no es suficiente para firmar archivos de métodos. También se precisa el derecho Editar métodos .			
Además, también se necesita el derecho Modificar métodos firmados para agregar la firma de aprobación a los archivos que ya tienen una firma de revisión.	Sí	No	No
Modificar métodos firmados*			
Editar un método firmado y guardarlo con el mismo nombre.	Sí	No	No
Ejecutar métodos no firmados*			
Con solo este derecho, únicamente se pueden usar los métodos aprobados.			
Nota: Este derecho se puede utilizar combinado con el derecho Ejecutar métodos no aprobados para permitir al usuario ejecutar todos los métodos.	Sí	Sí	No



	Administrador de magellan	Especialista en aplicaciones	Operador de magellan
Derechos del usuario		de magellan	g
Configurar favoritos			
Agregar/quitar metodos de la lista de favoritos.	Sí	No	No
Definir mediciones de etiquetas múltiples			
Definir métodos con mediciones de etiquetas múltiples.	Sí	Sí	No
Definir transformaciones cinéticas			
Definir métodos con transformaciones cinéticas.	Sí	Sí	No
Definir transformaciones de concentración			
Definir métodos con transformaciones de concentración.	Sí	Sí	No
Definir alias			
Definir métodos con seudónimos para los identificadores.	Sí	Sí	No
Zona de trabajo			
Crear zonas de trabajo			
Realizar una medición y guardar la zona de trabajo con un nuevo nombre, o bien, guardar una zona de trabajo existente con un nuevo			
nombre	Sí	Sí	Sí
Modificar zonas de trabajo			
Editar un espacio de trabajo ya existente y guardarlo con el mismo nombre.	Sí	Sí	No
Firmar zonas de trabajo*			
Firmar una zona de trabajo ya existente para revisión/aprobación.			
Nota: El derecho Firmar zonas de trabajo por sí solo no es suficiente para firmar archivos de zonas de trabajo. También se precisa el derecho Modificar zonas de trabajo .			
Además, también se necesita el derecho Modificar zonas de trabajo firmadas para agregar la firma de aprobación a los archivos que ya tienen una firma de revisión.	Sí	No	No
Editar zonas de trabajo firmadas*			
Editar un espacio de trabajo firmado y guardarlo con el mismo nombre.	Sí	No	No
Continuar evaluación en caso de error			
Continuar el cálculo después de producirse errores.	Sí	Sí	No ¹ /Sí ²

9. Icono Otras opciones



Derechos del usuario	Administrador de magellan	Especialista en aplicaciones	Operador de magellan
		de magenan	
Modificar datos sin procesar (muestras)			
muestras.	Sí	No	No
Modificar datos sin procesar (estándares)			
Editar o enmascarar datos de medición de			
estándares, controles, blancos.	Sí	No	No
Importar datos sin procesar			
Importar datos sin procesar de un archivo	Sí	No	No
Modificar evolucción		110	110
Cambiar los ajustes de evaluación (método)			
en una zona de trabajo existente.	Sí	Sí	No
Volver a evaluar con otro método			
Volver a evaluar una zona de trabajo			
existente mediante otro método.	Sí	Sí	No
Modificar diseño			
Cambiar los parámetros de medición, la concentración y el diseño antes de iniciar una			
medición	Sí	Sí	No
Curva estándar			
Crear archivos de curva estándar			
Cree una nueva curva estándar y guardarla con un nuevo nombre.	Sí	Sí	No
Editar archivos de curva estándar			
Editar una curva estándar existente y			
guardarla con el mismo nombre.	Si	Si	No
Firmar archivos de curva estándar*			
revisión/aprobación.	Reservado para uso futuro.		
Modificar archivos de curva estándar firmados*			
Editar y cambiar los ajustes de un archivo de			
curva estándar firmado.	Reservado para uso futuro.		
General			
Guardar en archivador			
Trasladar archivos a un archivador	Sí	No	No
Establecer directorios predeterminados**			
Establecer directorios individuales para			
guardar zonas de trabajo, metodos, listas de IDs de muestras, archivos exportados v			
mapas de bits.	Sí	Sí	Sí



Derechos del usuario	Administrador de magellan	Especialista en aplicaciones de magellan	Operador de magellan
Definir placas de filtro			
Cambiar la definición de filtros del instrumento.	Sí	No	No
Modificar opciones generales			
Elegir idioma, suprimir advertencia de no monótono, permitir modo demo.	Sí	No	No
Imprimir zonas de trabajo sin firmar*			
Se pueden imprimir los resultados de una zona de trabajo no firmada.	Sí	Sí	Sí
Ejecutar métodos no aprobados*			
Con solo este derecho se permite al usuario ejecutar métodos ya revisados, pero aún no aprobados, así como métodos aprobados. Nota: Este derecho se puede utilizar combinado con el derecho Ejecutar métodos no firmados para permitir al usuario ejecutar	Sí	Sí	No
todos los métodos.			
Editar método revisado* Los métodos revisados (pero no aprobados) se pueden editar.	Sí	No	No
Editar zona de trabajo revisada*			
Las zonas de trabajo revisadas (pero no aprobadas) se pueden editar.	Sí	No	No
Reader Server			
Guardar archivos de definición de placas			
Crear o editar archivos de definición de placas.	Sí	No ¹ /Sí ²	No
Guardar configuración de espín			
Crear o editar archivos de perfil de espín.	Sí	No	No

* derechos del usuario disponibles solamente en magellan Tracker

** derechos del usuario disponibles solamente en magellan Standard

¹ magellan Tracker

² magellan Standard



9.7 Acerca de magellan

Seleccione Acerca de magellan para ver la información de copyright y la exención de responsabilidad del software.

En el cuadro de diálogo **Acerca de magellan**, se muestra información acerca de la versión del software **magellan** instalada actualmente, los números de versión de los componentes (pestaña **componentes**), información UDI (Unique Device Identification) y la licencia del usuario del programa (pestaña **licencia**).

Asistente para registrar

Para conocer más detalles acerca del **Asistente para registrar magellan**, consulte el capítulo 2.4.4 Asistente para registrar.

10. Funciones adicionales de magellan Tracker

10.1 Administración de usuarios

En **magellan** Tracker, la administración de usuarios es obligatoria. Para obtener detalles adicionales, consulte el capítulo 9.4 Administración de usuarios (magellan Tracker).

10.1.1 Registro de modificaciones

Para cumplir las normas **FDA Regulation 21 CFR part 11** se documenta cada paso efectuado en **magellan** Tracker. En el **Registro de modificaciones** se enumeran todas las modificaciones relativas a los siguientes datos:

- un archivo de método
- un archivo de zona de trabajo
- una lista de IDs de muestras
- un archivo de curva estándar
- · la base de datos de administración de usuarios o
- opciones y ajustes predeterminados.

Cada anotación consta del usuario (nombre y nombre completo), fecha y la hora de la modificación, si el archivo ha sido creado o modificado, y los comentarios pertinentes.

Registro de modificaciones de Métodos, Zonas de trabajo y Listas de IDs de muestras

El Registro de modificaciones se puede consultar mediante el botón **Registro de modificaciones** de la página de guardar de los asistentes.

Registro de modificaciones de una Curva estándar

Las curvas estándar se pueden guardar y después se pueden utilizar para evaluar datos desde otra zona de trabajo. Para conocer los detalles de la función cargar una curva estándar desde un archivo externo, consulte la sección 4.3.8 Concentraciones: Curva estándar del Asistente Crear o editar método. Para abrir en el gráfico de curva estándar una curva estándar guardada previamente, consulte la sección 7.4.11 Barra de control: Concentraciones.

Para ver el registro de modificaciones de la curva estándar cargada, seleccione **Registro de modificaciones** en el menú **Ver** del cuadro de diálogo Gráfico de curva estándar.

Registro de modificaciones de la base de datos de administración de usuarios

Para ver el Registro de modificaciones de la base de datos de administración de usuarios, seleccione el icono **Otras opciones** en la página inicial de la **Lista de asistentes** y haga clic en **Administración de usuarios**.

Registro de modificaciones de Opciones y Configuración predeterminada

Para ver el Registro de modificaciones de todas las opciones y la configuración predeterminada, seleccione el icono **Otras opciones** en la página inicial de la **Lista de asistentes** y haga clic en **Opciones**. Entonces se podrá seleccionar en la pestaña **Otras opciones** el botón **Registro de modificaciones**.

10.2 Procesamiento de archivos

Los datos medidos siempre se guardan directamente después de la medición.

TECAN.



Aviso solo para Magellan Tracker: Para garantizar que no sea posible realizar ninguna manipulación (borrado) de los datos, los usuarios no deben ser administradores de Windows.

10.2.1 Guardar un archivo

Los archivos creados y guardados en **magellan** Tracker se señalan con un signo "#" en su nombre de archivo. El signo "#" añadido solamente se puede ver en el subdirectorio apropiado de **magellan**.

Ejemplo

método#1.mth

zonatrabajo#1.wsp

listalDmuestra#1.smp

10.2.2 Cambiar un archivo de zona de trabajo o método

Cada vez que se modifica y guarda un archivo de zona de trabajo o método, el número que sigue a "#" en el nombre de archivo se incrementa automáticamente (p. ej.,: primera versión de una zona de trabajo: nombre#1.wsp; edite y guarde la nueva versión para obtener nombre#2.wsp; Después del siguiente cambio se creará, nombre#3.wsp...). Todas las versiones de un archivo se enumeran en los subdirectorios correspondientes de **magellan**.



Aviso Cuando se transfieren archivos de Tracker es necesario transferir todas las versiones del archivo.

10.2.3 Abrir un archivo

Solamente la última versión guardada de un archivo está disponible en los asistentes de crear o editar un método, crear o editar una lista de IDs de muestras, evaluar resultados e iniciar medición.

Para ver una versión anterior de un archivo, abra el **Registro de modificaciones** en la página de guardar del archivo, haga clic una vez en el número de la versión que prefiera del archivo para seleccionarla y haga clic en el botón **Guardar versión previa como...** Cambie el nombre del archivo y añada algunas notas de archivo o comentarios de Registro de modificaciones y abra el archivo con el asistente **Crear o editar método**.



Aviso

Si desea comparar un método con sus versiones anteriores primero deberá imprimirlo, ya que no es posible abrir simultáneamente dos vistas preliminares de impresión.



10.2.4 Abrir un archivo creado en otro ordenador – Añadir HUIDs

En el proceso de asignación de licencias de **magellan**, hay tres números implicados:

- El número de serie del software
- El número de identificación de la unidad de hardware (HUID, Hardware Unit Identification Number) y
- El número de licencia (consulte el capítulo 2.4.3 Registrar magellan).

El HUID de un ordenador personal se configura durante el registro de **magellan** y todos los archivos creados y guardados en ese ordenador estarán vinculados con este HUID especial.

Si hubiera necesidad de trabajar con archivos de **magellan** Tracker creados en otro PC, se debe aceptar el HUID de este otro PC.

Cuando un usuario con derecho a **Modificar opciones generales** intenta abrir un documento (lista de IDs de muestras, método, zona de trabajo) desde otro PC, es decir, con un número HUID distinto, al usuario se le pedirá automáticamente que añada ese número a la lista de números HUID conocidos. Si se hace clic en el botón **Aceptar HUID**, ese número HUID se añadirá a la lista y se abrirá el documento sin interacciones adicionales.

En lo sucesivo ese número HUID siempre será aceptado.

Para ver la lista de números HUID aceptados de otros PCs, haga clic en el icono **Otras opciones** en la página de inicio de la lista de asistentes \rightarrow **Opciones** \rightarrow pestaña **Otras opciones**. Haga clic en el botón **Añadir HUID** y se mostrará el cuadro de diálogo que contiene los números HUID de otros PCs.

HUID no válido	×
Este archivo fue creado en otro ordenador (HUID: 9408G-3U335FB). (C: \Users\Public\Documents\Tecan\Magellan\mth\Application Example_en	n.mth)
Pulse 'Aceptar HUID' para agregar el HUID dado a la lista de HUIDs dispo lista se pueden abrir sin recibir ningún mensaje de advertencia.	nibles. Los archivos de los ordenadores de esta
	Aceptar HUID Rechazar HUID

10.3 Registro de modificaciones del sistema

El registro de modificaciones del sistema solo está disponible en la versión **magellan Tracker**. La función de registro de modificaciones del sistema escribe varios eventos de **magellan** en un archivo de registro (véase la siguiente tabla). Una vez por semana se crea un nuevo archivo de registro y se guarda en

Windows 10: C:\Users\Public\Documents\Tecan\LogFiles\SystemAuditTrail

Se usa una suma de control calculada sobre el archivo de registro completo para garantizar la consistencia del archivo de registro y proteger los archivos frente a cambios fortuitos e intencionados. Durante el inicio siempre se comprueba la coherencia del archivo de registro. Los archivos no válidos se cerrarán inmediatamente, se marcarán como no válidos anexando ".invalid" al nombre del archivo de registro y se creará uno nuevo en su lugar.

Los archivos de registro antiguos o no válidos se mueven a la carpeta de Archivo, donde están protegidos contra la eliminación.

Evento	Tipo de evento	Descripción
Iniciar magellan	Iniciar	Se inició magellan.
Salir de magellan	Salir	Se cerró magellan.
Inicio de sesión de usuario correcto	Inicio de sesión	Un usuario ha iniciado sesión correctamente.
Inicio de sesión de usuario fallido	Inicio de sesión erróneo	Inicio de sesión fallido porque el nombre de usuario o la contraseña son incorrectos.
Cierre de sesión de usuario	Cerrar sesión	Un usuario ha cerrado sesión.
Nueva zona de trabajo	Nueva	Se creó una nueva zona de trabajo.
Cargar zona de trabajo	Cargar	Se cargó en magellan un archivo de zona de trabajo existente.
Guardar área de trabajo	Guardar	Se guardó una zona de trabajo.
Nuevo método	Nuevo	Se creó un nuevo método
Cargar método	Cargar	Se cargó en magellan un archivo de método existente.
Guardar método	Guardar	Se guardó un método.
Nueva lista de IDs de muestras	Nueva	Se creó una nueva lista de IDs de muestras.
Cargar lista de IDs de muestras	Cargar	Se cargó una nueva lista de IDs de muestras.
Guardar lista de IDs de muestras	Guardar	Se guardó una lista de IDs de muestras.
Iniciar medición	Medir	Se inició una nueva medición.
Adjuntar firma	Firmar	Se firmó un archivo de zona de trabajo o de método.
Exportar resultados	ExportASCII ExportExcel ExportASTM ExportSampleTra cking	Se exportaron los resultados.
Imprimir resultados	Imprimir	Se imprimieron los resultados.
Archivado automático	ArchiveFailed	Error al archivar un archivo de zona de trabajo, método o lista de IDs de muestras.
Error de CRC	Fallo de CRC	La suma de control del archivo de registro anterior no era válida.

Eventos de magellan de los que se hará seguimiento

11. Cálculos

11.1 Evaluar resultados – Procedimiento de cálculo

En función de la configuración, el cálculo sigue este procedimiento:

- 1. Cálculo previo
 - 1.1 Cálculo de espectros
- 2. Comprobar si encajan los datos y la configuración de la evaluación
- 3. Cálculo de estadísticas de datos sin procesar
- 4. Cálculo de transformación
- 5. Cálculo de parámetro cinético
- 6. Cálculo de transformación cinética
- 7. Cálculo de concentración
- 8. Cálculo de transformación de concentración
- 9. Determinación del rango de corte

10.Validación QC

11.Cálculo estadístico

El procedimiento se repite si

- La validación de QC eliminó valores mediante funciones de eliminación.
- · Las transformaciones requieren parámetros de interpolación.

Se suprimirán los valores iguales para transformaciones en cinéticas y en mediciones de etiquetas múltiples.

Para los cálculos con diluciones diferentes, se desecharán las concentraciones Medias.

En caso de que se interrumpa el procedimiento de evaluación debido a errores, se mostrarán los resultados calculados hasta ese punto.

En caso de que algunos valores resulten ser FALSOS, se mostrará el siguiente cuadro de diálogo **Detectados errores**:

Campo de texto del mensaje	En un campo de texto se muestran el mensaje de error y una breve explicación.
Botón Guardar como archivo	Haga clic en el botón Guardar como archivo y podrá guardar en un archivo ASCII el protocolo de error.

En caso de que falle el ajuste de curva, dependiendo de los derechos del usuario, se mostrará una pregunta para que el usuario pueda decidir cambiar los ajustes del Tipo de análisis. Después de hacer los cambios, el procedimiento de evaluación se iniciará desde el principio.

En caso de que falle la validación, dependiendo de los derechos del usuario, se mostrará un cuadro de mensaje de error o un cuadro de mensaje de pregunta, donde el usuario podrá decidir si desea continuar el cálculo de todos modos. Si se interrumpe el cálculo, no se podrá tener acceso a los resultados de corte. No obstante, en la copia impresa se incluirá un mensaje de fallo de validación.

En caso de que se hayan cambiado los valores o los ajustes, se mostrará un **cuadro de mensaje**.

Mensaje	Han cambiado los valores del instrumento. ¿Desea reiniciar el cálculo resultante?
Botón Sí	Haga clic en el botón S í para calcular de nuevo los valores de evaluación.
Botón No	Haga clic en el botón No para cerrar ese mensaje. No se realizará ningún recálculo.

11.1.1 Estadísticas

Si se calcula un valor estadístico p. ej., sobre las n réplicas con valores $x_1...x_n$ de una muestra, se utilizan las siguientes fórmulas:

Nombre	Fórmula
Valor medio, valor promedio	$\overline{x} = \frac{1}{n} \sum_{i=1}^{n} x_i$
Desviación estándar	$s = \sqrt{\frac{1}{n-1}\sum_{i=1}^{n} (x_i - \bar{x})^2}$ y s=0 para n=1
Coeficiente de variación, en porcentaje	$v = 100 \frac{s}{\overline{x}}$

Para obtener más información, consulte la sección 11.3.5 Funciones estadísticas.

11.2 Reducción de datos de espectros

Los datos de espectros contienen mucha más información que simples mediciones de longitudes de onda. Entre la información que habitualmente se extrae de los espectros están las intensidades a longitudes de onda especificadas o en picos encontrados. Durante el cálculo también se pueden calcular la reducción del blanco espectral y diversas relaciones.

11.2.1 Descripción matemática

La **reducción del blanco** se lleva a cabo restando el espectro blanco del espectro de muestra.

El **suavizado** se realiza con el algoritmo de Savitzky-Golay. Los coeficientes del filtro se calculan con un polinomio de grado 2. El factor de suavizado se trata como número de puntos:

$$g_i = \sum_{n=-n_i}^{n_R} c_n f_{i+n}$$
 donde n_R-n_L+1=número de puntos

La ventana de suavizado es simétrica, p. ej., para un factor de alisamiento de 5, n_L =-2, n_R =2.

Para los bordes inferior y superior del espectro, la ventana es asimétrica, así pues, $n_L=0$, $n_R=4$ o $n_L=-4$, $n_R=0$ respectivamente.

Antes del suavizado, se eliminan del espectro todos los valores de desbordamiento.

Selección de longitud de onda





- Seleccionar longitud de onda Obtiene la intensidad de la longitud de onda seleccionada
- Calcular proporción Calcula la relación dividiendo las intensidades a las longitudes de onda seleccionadas
- Área Calcula el área bajo el espectro dentro del rango seleccionado de longitudes de onda
- A definir por el usuario Reduce el espectro utilizando una fórmula definida por el usuario

Normalizar

Reduce el espectro en un rango de 0 a 1.

Encontrar pico

Se diferencia el espectro y todos los puntos nulos del resultado se tratan como picos potenciales. Los picos correspondientes se ordenan por valor y todos los picos no válidos se eliminan (p. ej., desbordamientos, mínimos locales, valores por debajo del umbral definido). El valor más alto de los elementos restantes se selecciona como pico.

En la práctica, este es a menudo el valor máximo del espectro.

Fórmulas personalizadas

Capacidad de definir diversas reducciones de datos para los espectros (consulte la sección 11.3.8 Funciones de espectros).



ADVERTENCIA

LOS MARGINALES AISLADOS DEBIDOS AL RUIDO TAMBIÉN SE DETECTARÁN COMO PICOS. ¡POR TANTO SE RECOMIENDA ENCARECIDAMENTE SUAVIZAR EL ESPECTRO ANTES DE UTILIZAR ENCONTRAR PICO!

11.3 Escribir una fórmula

11.3.1 Introducción

Los cálculos, los requisitos de corte y la validación requieren diversas fórmulas. Estas fórmulas tienen una sintaxis semejante a los comandos de BASIC.

Todos los valores se calculan con doble precisión, aunque se muestren en el formato numérico deseado.

11.3.2 Variables en fórmulas

Todas las abreviaturas definidas previamente para los identificadores contenidos en las placas de análisis se pueden emplear como variables en los cálculos.

Por ejemplo:

PC1, SM1_1, BL1...Valores promedio

PC1_1, SM1_1_1, BL1_1...Valores individuales, réplicas individuales



Aviso Cuando procesa las etiquetas, el programa diferencia entre letras mayúsculas y minúsculas.

También se puede trabajar con coordenadas de pocillo.

Por ejemplo: A2, B3, H12





Aviso El símbolo x se refiere al valor actual dentro de un pocillo. Por ejemplo: x-BL1...resta el valor vacío (blanco) del valor actual.

Aviso Las coordenadas de pocillo siempre se deben escribir con letras mayúsculas.



Aviso Se puede tener acceso a los distintos ciclos para mediciones cinéticas por medio de índices [], donde"0" indica el primer ciclo.

Por ejemplo:

SM1_1[0]... Valor promedio de la muestra 1 del primer ciclo

x[1]... valor de pocillo actual del segundo ciclo

x[i] ... valor del pocillo actual y el ciclo actual

Substracción cinética consecutiva – Cálculo de diferencias entre ciclos cinéticos

Se pueden definir cálculos sobre datos cinéticos mediante la variable de iteración 'i' en las fórmulas. Por ejemplo, se puede realizar una derivada numérica de una curva cinética mediante la fórmula x[i]-x[i-1], que resta cada ciclo cinético del anterior. Pero el uso de esa fórmula conduciría automáticamente a un error para el ciclo número 1. Por lo tanto se debe usar la fórmula **if(i>0) then(x[i]-x[i-1]) else ignore()** en el campo de edición de la transformación para evitar la substracción del primer ciclo.



El uso de constantes en las fórmulas es idéntico a todas las demás variables, la única diferencia es que una constante puede contener caracteres (p. ej., alfa, dilución).

Hay dos variables de acceso predefinidas.

concX ...

devuelve la concentración del estándar del pocillo actual

dilX …

devuelve la dilución de la muestra o el control del pocillo actual

Aviso



Si se va a utilizar más de un conjunto de datos de entrada para los cálculos, seleccione los datos apropiados en la lista desplegable de datos disponibles, activados en la pestaña de los campos de editar transformación de Crear o editar método (consulte el capítulo 4.3.5 Datos transformados: Añadir nueva transformación y el capítulo 7.4.9 Barra de control: Datos transformados).



Aviso Al utilizar las funciones matemáticas que se describen en las secciones siguientes, se activará una ayuda de tipo "IntelliSense" para apoyar la escritura correcta de fórmulas especiales.

Operadores de fórmulas

Operador	Descripción
+	Más
-	Menos
*	Multiplicación
1	División
^	Elevado a la potencia del exponente
	El símbolo utilizado para representar la expresión "a la potencia de" es '^'. En un teclado inglés, se puede pulsar la tecla Mayúsculas junto con la tecla número 6 (no en el teclado numérico) para obtener ese carácter. Ejemplo: x^3 El valor del pocillo se elevará a la tercera potencia.
<	Menor que
<=	Menor o igual que
>	Mayor que
>=	Mayor o igual que
==	Igual a
!=	No igual a
=	Asignado a





Aviso Hay diferencias entre igual a (==) y asignado a (=). Por ejemplo: If (x == 0,000), then x = 1,000 (Si el valor actual de x es igual a 0 entonces asignar a x el valor 1)

11.3.3 Funciones de fórmulas

Las expresiones "and" y "or" se pueden incorporar a la misma ecuación a través del uso de funciones lógicas. El resultado de una ecuación lógica será VERDADERO o FALSO y por lo tanto es un valor booleano. Las ecuaciones individuales se deben encerrar entre paréntesis.

Expresiones lógicas

and

En ecuaciones de esta naturaleza, el resultado será VERDADERO si ambas expresiones resultan ciertas; en caso contrario, darán como resultado FALSO. Ejemplo:

Si el valor que tiene el pocillo es 0,3

(x>0,0) and (x<1) = TRUE (x>0,0) and (x<0,1) = FALSE

or

En ecuaciones de esta naturaleza, se debe cumplir que al menos una de las expresiones implicadas es cierta para que se produzca el resultado VERDADERO.

Ejemplo: si el valor que tiene el pocillo es 0,3

(x>0,0) or (x<0,1) = TRUE (x>0,4) or (x<0,1) = FALSE

if(...) then(...) else(...)

Las siguientes expresiones se utilizan en el programa y, como ejemplo, se pueden escribir de la siguiente manera:

if (expresión) then Término A else Término B

La expresión:

la expresión debe ser verdadera o falsa y se debe escribir entre paréntesis. Si la expresión se obtiene por medio de operadores lógicos (and/or), entonces toda la expresión se debe escribir aquí y se debe encerrar entre paréntesis.

Términos A y B:

si la expresión es verdadera, se obtendrá el Término A y si la expresión es falsa se obtendrá el Término B.

No siempre es necesario incluir el operador else. Si no está presente en la ecuación, no se dará ninguna respuesta en caso de un resultado falso.

La expresión "if(...) then(...) else(...)" también se puede utilizar con términos booleanos para la validación QC.
Ejemplo:

Comprobar si el valor de un pocillo está comprendido dentro de un rango definido. Si el resultado es sí, el pocillo se representará por el valor 0,0 y si el resultado es no, se usará el valor 1.

if ((x>-0,005) and (x<0,0)) then(0,0) else(1,0)

11.3.4 Funciones básicas

abs(argumento)	
	Esta función devuelve la suma absoluta del argumento.
	Es decir, devuelve (-1*x) cuando x<0 y (x) cuando x>=0.
	Ejemplo:
	abs(-1) = 1
	abs(1) = 1
exp(argumento)	
	Esta función eleva el número de Euler (e) a la potencia del argumento.
	Ejemplo: $exp(1) = e^1 = 2,718$
frac(argumento)	
	Esta función separa la parte fraccionaria del argumento.
	frac(): devuelve la parte fraccionaria de un valor decimal.
	Ejemplo: $frac(1,7) = 0,7$
int(argumento)	
	Esta función separa la parte entera del argumento
	int(): devuelve la parte entera de un valor decimal
	Eiemplo: <i>int</i> $(1,7) = 1$
Log(argumento)	log(arg1:arg2)
	Esta función toma el logaritmo de arg2 signdo la base arg1
	Eigendo:
	El posillo tiono un volor de 100
	$\log(x, TO)$
	El resultado de este ejemplo sera 0,5.
In(argumento)	
	Esta función calcula el logaritmo natural del argumento.
	Ejemplo : <i>In</i> (10) = 2,303
lg(argumento)	
	Esta función calcula el logaritmo en base 10 del argumento.
	Ejemplo: $lg(10) = 1$



round(argumento)

Esta función redondea el argumento a un valor entero.

Ejemplo:

Para redondear 12,579 a dos decimales se puede utilizar la siguiente fórmula:

round(12,579*100)/100

El resultado es 12,58.

sqr(argumento)

Esta función eleva el argumento al cuadrado.

Ejemplo: sqr(3) = 9

sqrt(argumento)

Esta función calcula la raíz cuadrada del argumento.

Ejemplo: sqrt(9) = 3

11.3.5 Funciones estadísticas



Aviso Para transformaciones con funciones estadísticas, utilice datos simples como datos de entrada, no promedios ni medias.

avg(argumento) y mean(argumento)

Esta función calcula el valor promedio del argumento. El argumento debe ser un identificador.

Ejemplo:

avg(SM1)

Se obtendrá el promedio de todas las muestras del primer grupo experimental. Además, se puede calcular el promedio de las réplicas de un valor:

avg(ST1_1)

Promedia todas las réplicas del primer estándar.

median(argumento)

Esta función calcula la mediana del argumento. El argumento debe ser un identificador. Las réplicas individuales se ordenan según su tamaño y la mediana se determina tomando el valor medio en esa ordenación. Si el número de valores es un número par, se determinará el promedio de los dos valores medios.

Ejemplo:

median (NC1)
Supongamos que NC1_1=0,1, NC1_2=0,05, NC1_3=0,04
El valor de la mediana de estos controles negativos se determinará como sigue.
Al ordenar los valores, devuelve:
0,04 0,05 0,1
La mediana es por lo tanto 0,05.



medianPlate()

Esta función se utiliza para determinar la mediana de la placa completa. Los valores individuales se ordenan según su tamaño y la mediana se determina tomando el valor medio en esa ordenación. Si el número de valores es un número par, se determinará el promedio de los dos valores medios.

Ejemplo:

medianPlate()

Supongamos que BL1=1, NC1=2, PC1=3, SM1_1=4, SM1_2=5 y que no hay definido ningún otro pocillo.

La mediana de la placa completa es 3.

max(argumento) y min(argumento)

Esta función calcula el máximo/mínimo del argumento. El argumento debe ser un identificador.

Ejemplo:

max(SM1)

Se obtendrá el máximo de todas las muestras del primer grupo experimental.

min(ST1_1)

Determina mínimo de las réplicas individuales del primer estándar.

maxAvg(argumento) y minAvg(argumento)

Esta función devuelve el valor promedio máximo/mínimo del argumento. El argumento debe ser un identificador.

Ejemplo:

minAvg(SM1)

Si para SM1 hubiera varios resultados, solamente se devolverá el promedio mínimo.

PointwiseCV(argumento)

Determina el coeficiente de variación (CV) promedio para cada punto. El resultado se puede utilizar para validar la curva estándar. El argumento debe ser un identificador.

Ejemplo:

PointwiseCV(ST1)

Supongamos que

ST1_1_1=0,54 ST1_1_2=0,52 cv=2,668

ST1_2_1=0,72 ST1_2_2=0,77 cv=4,746

ST1_3_1=1,08 ST1_3_2=0,99 cv=6,148

El resultado de pointwiseCV es el promedio del CV de ST1_1, ST1_2 y ST1_3, que resulta 4,5209.

Sum(argumento)

Esta función calcula la suma del argumento.

Ejemplo:

sum(SM1)

Se obtendrá la suma de todas las muestras del primer grupo experimental.

sum(ST1_1)

Determina la suma de las réplicas del primer estándar.

stddev(argumento)

Esta función calcula la desviación estándar del argumento. El argumento debe ser un identificador.

Ejemplo:

stddev(NC1)

Se obtendrá la desviación estándar del control negativo del primer grupo experimental.

11.3.6 Funciones de eliminación

Las funciones de eliminación se emplearán en la validación. De esta manera el usuario puede estar seguro de que los valores de medición están comprendidos en un intervalo válido y eliminar cualquier valor inapropiado.

Al definir los intervalos de validez, se pueden emplear tres funciones de eliminación distintas. Cuando se definen los criterios de eliminación, el primer parámetro requerido siempre es el identificador al que va a afectar la función.

Las tres funciones de eliminación producen un resultado lógico. Devolverá VERDADERO cuando haya suficientes valores de medición válidos disponibles y FALSO cuando haya menos que los requeridos.

Los valores de medición no válidos serán etiquetarán con un signo de admiración.

eliminate (arg1; arg2; arg3; arg4)

Se definirá el intervalo de validación mediante un valor definido.

Esta función elimina los valores que están fuera del rango de validación definido en arg2.

arg1: nombre de identificador (NC1, PC1).

arg2: valor utilizado para representar el rango, depende de arg4.

arg3: número de valores válidos individuales que deben estar disponibles para poder generar un promedio. Si quedan pocos valores disponibles después de la eliminación, se devolverá el resultado de FALSO.



Aviso Para arg3: se debe definir un valor de "2" o superior. "1" no es aceptable.

arg4: representa la selección de si la prueba utilizará el rango promedioarg2 a promedio+arg2 (argumento 1) o si se comprobará si los valores individuales se encuentran por debajo de promedio+arg2 (argumento 0). En este caso (argumento 0) no hay ningún límite inferior.

La función de eliminación aquí determina el promedio de arg1. A continuación se compara el valor con la mayor desviación absoluta respecto al valor promedio



para asegurarse de que está comprendido en el rango que indica arg4. Si el valor se encuentra fuera de ese rango, se considerará no válido. En ese caso, el cálculo del promedio comenzará de nuevo excluyendo el valor que previamente se descubrió que era no válido y se continuará de este modo hasta que no se pueda encontrar ningún valor inapropiado más.

Se devolverá un resultado de VERDADERO cuando haya disponibles suficientes valores individuales válidos.

Ejemplo:

eliminate(NC1;0,15;2;0)

Se generará el valor promedio para los controles negativos. Se comprobará a continuación si los controles negativos individuales se encuentran por encima de NC1+0,15. (No hay ningún límite inferior). Si es así, se eliminarán los valores inapropiados como se ha descrito. Si al finalizar, al menos dos valores individuales son válidos, se devolverá el resultado VERDADERO; de lo contrario, se devolverá el resultado FALSO.

eliminatePerc (arg1; arg2; arg3; arg4)

Se definirá el intervalo de validación mediante un porcentaje del valor promedio.

Esta función elimina valores como se ha descrito para la función eliminate. El rango de validación se define mediante un porcentaje (arg2) del identificador.

Ejemplo:

eliminatePerc(NC1;10;2;1)

Los valores individuales serán no válidos si se desvían hasta una amplitud de más del 10% por debajo o por encima del promedio. En cuanto se termine, deben ser válidos un mínimo de dos valores individuales para que los controles negativos se declaren válidos.

eliminateRange (arg1; arg2; arg3; arg4)

Esta función elimina los valores comprendidos fuera del rango de validación definido por los argumentos 2 y 3.

arg1: nombre de identificador (NC1, PC1)

arg2: Límite inferior del rango seleccionado

arg3: Límite superior del rango seleccionado

arg4: Es el número de valores individuales válidos necesarios para continuar. Si quedan pocos valores disponibles después de la eliminación, se devolverá el resultado de FALSO.

Se devolverá un resultado de VERDADERO cuando haya disponibles suficientes valores individuales válidos.

Ejemplo:

eliminateRange(NC1; 0,0; 0,1; 2)

Para que sean válidos, los valores individuales para los controles negativos deben estar comprendidos dentro del rango de 0,0 a 0,1. Para conseguir un resultado VERDADERO, al menos dos de los valores se deben considerar válidos.



eliminateCV(arg1; arg2; arg3)

Esta función elimina réplicas hasta que el CV de las réplicas restantes sea inferior que el CV dado (arg2). Las réplicas se eliminan paso a paso, comenzando por el valor con la mayor diferencia respecto al valor medio.

arg1: Nombre de identificador

arg2: Límite del valor de CV aceptado

arg3: Ese el número de valores individuales válidos que deben quedar para que se devuelva un resultado positivo. Si quedan pocos valores disponibles después de la eliminación, se devolverá el resultado de FALSO.

Se devolverá un resultado de VERDADERO cuando haya disponibles suficientes valores individuales válidos.

Ejemplo:

eliminateCV(NC1; 15; 3)

El CV calculado de los valores individuales del control negativo deberá estar por debajo del 15 %. Después de la eliminación, deben quedar al menos tres réplicas para obtener VERDADERO como resultado.

countDeleted(arg1; arg2)

Esta función comprueba si hay suficientes valores válidos disponibles y devuelve VERDADERO o FALSO.

arg1: El identificador, cuya cantidad se debe comprobar.

arg2: El menor número de réplicas que deben estar disponibles.

Ejemplo:

countDeleted(NC1; 2)

Esta función trabaja en las ecuaciones junto con las funciones de Eliminar:

La Validación contiene la siguiente línea:

if (NC1>0,5) then eliminateRange(NC1; 0; 0,5; 2)

Si el promedio del control negativo se encuentra por encima de 0,5, entonces se deberán omitir todas las réplicas comprendidas fuera del rango.

Esta eliminación se procesará una vez. Si se encuentran valores para eliminar, seguirá un nuevo cálculo para asegurarse de que el promedio se encuentra por debajo de 0,5 y se devolverá un resultado de VERDADERO.

En ese punto todavía no se está seguro de si está disponible el número necesario de réplicas. Entonces se determinará ese detalle mediante la función countDeleted dentro de una expresión else:

if (NC1>0,5) then eliminateRange(NC1; 0; 0,5; 2) else countDeleted(NC1; 2)

El proceso de ecuación (de la segunda ejecución después de la eliminación) utiliza la expresión else. La función countDeleted entonces comprueba, si el identificador devuelto en arg1 está disponible en la cantidad especificada en arg2. Si es así, se devolverá un resultado de VERDADERO, o FALSO en caso contrario.

Ejemplo paso a paso del uso de eliminate (también válido para eliminatePerc)

La única diferencia entre eliminate y eliminatePerc es el modo de calcular los límites para el rango válido (consulte los párrafos anteriores). La secuencia de cálculo es la misma para ambas funciones. En la descripción paso a paso, se utiliza eliminate.

eliminate(NC1;0,15;2;0)

Valores individuales:	NC1_1 = 0,217	
	NC1_2 = 0,439	$\overline{NC1} = 0.288$
	NC1_3 = 0,208	

Paso 1: se ordenan los valores individuales según su distancia al promedio.

$ NC1_1 - \overline{NC1} $	0.217-0.288	0.071		NC1_2 (0,439)
$ NC1_2 - \overline{NC1} $	0.439-0.288	0.151	\rightarrow	NC1_3 (0,208)
$ NC1_3 - \overline{NC1} $	0.208-0.288	0.08		NC1_1 (0,217)

Paso 2: cálculo del límite superior, 0.288+0.15 = 0.438

Paso 3: primera comparación:

0,439 > 0,438 y por lo tanto fuera del rango válido \rightarrow Paso 4: se elimina el valor de NC1_2 y se calcula de nuevo el promedio y el límite.

Paso 4: Nuevo promedio = (0,217+0,208)/2 = 0,2125

Nuevo límite = 0,2125+0,15 = 0,3625

Paso 5: segunda comparación:

0.208 <= 0.3625 → válido

0.217 <= 0.3625 → válido

Resultado:

La réplica NC1_2 no encaja en el rango devuelto y por lo tanto ha sido eliminada.

La función devuelve VERDADERO porque todavía quedan dos réplicas.

Aviso: En cada ciclo de cálculo (Paso 1 a Paso 4), solamente se elimina una réplica (la réplica con mayor distancia al promedio). El ciclo se repite hasta que no se elimine ninguna réplica durante el paso de comparación dentro del ciclo.

11.3.7 Otras funciones

ignore()

Esta función omite los valores de los pocillos seleccionados para no sean utilizados en los cálculos. Los valores que no se tienen en cuenta se muestran precedidos por el signo '!' y se procesan como valores eliminados.

Ejemplo:

if(x<0,9*SM1_1) then ignore()

isInvalid()

Esta función devuelve un valor booleano. Indica si los valores dentro del pocillo son válidos o no. Los valores no válidos, por ejemplo, pueden estar ocasionados por valores de desbordamiento del lector.

Ejemplo:

if(isInvalid()) then x=3,0

Si el pocillo contiene un valor no válido, se le asignará el valor 3,0. De esta manera, será posible realizar cálculos adicionales, a pesar de tales valores.

calcAlways(argumento)

Esta función permite calcular fórmulas sobre pocillos independientemente del estado del valor. El argumento puede ser cualquier fórmula válida. La fórmula se calculará aunque el valor esté enmascarado o sea no válido. Esta función se puede utilizar para calcular resultados que no dependen del valor del pocillo utilizado.

Ejemplo:

calcAlways(A)

Devuelve el valor del parámetro A de la fórmula de curva estándar.

concX

Esta función solamente se puede utilizar en transformaciones de concentración. El resultado de **concX** es la concentración original del pocillo y por lo tanto solamente se puede utilizar en pocillos con estándares como identificador.

Ejemplo:

x – concX devuelve la diferencia entre la concentración calculada y original del pocillo.

11.3.8 Funciones de espectros

smooth(númPts)

Suaviza con el número dado de puntos.

minimum()

Devuelve el mínimo como un par de datos (longitud de onda, intensidad).

maximum()

Devuelve el máximo como un par de datos (longitud de onda, intensidad).

normalize(lo)

Normaliza respecto a la intensidad para la longitud de onda dada.

peak(umbral;loInicial;loFinal;índicePico:porcLímite)

Devuelve el pico como un conjunto de datos (intensidad, longitud de onda, anchura, área). No se encontrarán los picos inferiores al umbral. La búsqueda del pico se realiza en el rango dado de longitudes de onda.

índicePico = $-1 \rightarrow$ devuelve el pico máximo

índicePico = $-2 \rightarrow$ devuelve el pico mínimo

índicePico = 0, 1, 2, ... \rightarrow devuelve el pico para el índice dado

porcLímite es un porcentaje utilizado para calcular la anchura y el área.

numPeaks(umbral;loInicial;loFinal)

Devuelve el número de picos encontrados en el rango dado de longitudes de onda. No se encontrarán los picos inferiores al umbral.

derive(grado)

Calcula la deriva del grado dado. Los grados válidos son 1 y 2; p. ej., derive(2).

intensity(función)

Devuelve la intensidad de un par de datos o de un pico. P. ej., intensity(minimum()).



wavelength(función)

Devuelve la longitud de onda de un par de datos o de un pico. P. ej., wavelength(maximum())

width(función)

Devuelve la anchura de un pico; p. ej., width(peak(0,5;450;650;-1;50)).

area(función)

Devuelve el área de un pico; p. ej., area(peak(0,5;450;650;-1;50)).

11.3.9 Ejemplos

Transformaciones

Las transformaciones modifican el valor actual del pocillo. El resultado siempre debe ser un valor numérico.

Reducción del blanco:

x-BL1

Se deducirá el valor vacío (blanco) del valor actual del pocillo.

Blanking de una ejecución cinética:

x-x[0]

El valor del primer punto de la cinética se utiliza como blanco y se deducirá de los valores de todos los ciclos cinéticos.



Aviso

El índice "0" representa la primera medición de una ejecución cinética.

Proporción:

x/PC1 * 100

Calcula el cociente entre una muestra y un control en porcentaje.

Proporción ADN/proteínas:

'Etiqueta1'!x/'Etiqueta2'!x

Para calcular la proporción de ADN/proteínas, se debe definir una medición de etiquetas múltiples. La primera medición utiliza una longitud de onda de 260 nm y la segunda medición utiliza una longitud de onda de 280 nm.

Resultado condicional:

If (x<0,0) then 0 else x

Si el valor del pocillo es menor que 0, será representado por 0, de lo contrario permanecerá sin cambios.

Corte

El corte dicta los límites que se aplicarán al evaluar. El resultado siempre debe ser un valor numérico.

Etiqueta: POS Límite: NC1*1,15 Etiqueta: ?? Límite: NC1*0,85 Etiqueta: -



El resultado será positivo (POS) cuando la absorbencia esté más del 15% por encima del promedio de los controles negativos, y será negativo (-) cuando esté más del 15% por debajo del promedio. Un resultado se declara abierto (??) si se encuentra entre esos dos valores.

Validación QC

Esta validación sirve para examinar la validez de la prueba. El resultado siempre debe ser una expresión lógica, VERDADERO o FALSO.

(BL1>-0,005) and (BL1<0,120)

Si el valor promedio del valor vacío está comprendido dentro del rango -0,005 a 0,12, se declarará un resultado de VERDADERO.

eliminateRange(NC1;-0,005;0,12;2)

Si un mínimo de dos controles negativos están comprendidos dentro del rango - 0,005 a 0,12, se declarará un resultado de VERDADERO.

abs(PC1-NC1)>0,25

La distancia entre los promedios de los controles positivos y los controles negativos debe ser mayor que 0,25. Si es así, se declarará un resultado de VERDADERO.

11.4 Tipos de análisis de la curva estándar

11.4.1 Definiciones

Datos n puntos base $(x_1, y_1), ..., (x_n, y_n), x_1 < ... < x_n$

Nombre	Fórmula
Coeficiente de correlación	$r = \frac{\sum_{i=1}^{n} (x_i - \bar{x})(y_i - \bar{y})}{\sqrt{\sum_{i=1}^{n} (x_i - \bar{x})^2 \sum_{i=1}^{n} (y_i - \bar{y})^2}}$
Desviación cuadrática media	$d = \sqrt{\frac{1}{n} \sum_{i=1}^{n} (y_i - f(x_i))^2}$
Calidad del ajuste	goodness = 1 - d

11.4.2 Parámetros de tipo de análisis

En el cálculo se puede usar el parámetro de tipo de análisis (p. ej., pendiente A e intercepción B para la regresión lineal). A continuación se describen los parámetros disponibles. **Si hay más de un grupo experimental**, los parámetros de tipo de análisis tienen un sufijo "B" para el segundo grupo experimental, "C" para el tercer grupo experimental, etcétera. Ejemplo: a la pendiente de una regresión lineal en el grupo experimental 3 se puede acceder como "AC".



11.4.3 Mensajes de error

Si cualquiera de los valores en las concentraciones calculadas se encuentra fuera del rango especificado por los estándares, se mostrará, respectivamente, la entrada >Máx. o <Mín. dentro de los pocillos de la placa. No obstante, si el usuario quisiera incorporar estos valores al cálculo, debe seleccionarse la opción Extrapolación.

Si la curva no es estrictamente monótona, se generará un mensaje de error.

Si hay varias soluciones para un cálculo de concentración para un valor medido, se mostrará la entrada MultPt dentro de los pocillos de la placa.

11.4.4 Punto a punto

Utilizando este proceso, los puntos base adyacentes se unirán por medio de una línea recta.

Este cálculo requiere un mínimo de 2 puntos base.

La extrapolación no es posible. En las transformaciones no hay ningún parámetro que pueda usarse como variable.

Si no se puede calcular ninguna concentración, se mostrará la entrada NoCalc dentro de los pocillos de la placa.

Descripción matemática

Función de interpolación:

$$f:[x_1, x_n] \to R$$

$$x \mapsto y_i + \frac{y_{i+1} - y_i}{x_{i+1} - x_i} \cdot (x - x_i) \quad \text{if } x \in [x_i, x_{i+1}]$$

11.4.5 Regresión lineal

Se formará una línea recta donde la suma de las desviaciones cuadráticas de todos los puntos bases es un mínimo.

Este cálculo requiere un mínimo de 2 puntos base.

Si fuera necesario, también se puede implementar extrapolación en este proceso. En una evaluación con un grupo experimental, la pendiente y la intercepción de ordenada sirven como variables A y B en las transformaciones. Además, también se puede implementar el coeficiente de correlación r.

En el cuadro de diálogo Curva estándar se muestra la información adicional siguiente:

Fórmula con valores de coeficientes A (pendiente) y B (intercepción)

Desviación cuadrática media d

Valor absoluto del coeficiente de correlación r

Descripción matemática

Función de aproximación:

$$f:[x_1,x_n]\to R$$

$$x \mapsto A \cdot x + B$$

donde A y B se determinan minimizando la función de error

$$err(A, B) = \sum_{i=1}^{n} (f(x_i) - y_i)^2$$

La solución es única si

$$rank \begin{pmatrix} x_1 & 1 \\ \vdots & \vdots \\ x_n & 1 \end{pmatrix} = n$$

que es verdadero si $x_i \neq x_j \forall i, j = 1,...,n$ (véase condición general)

11.4.6 Regresión no lineal

Esta aproximación está diseñada especialmente para datos hiperbólicos. Idealmente, los valores medidos para concentraciones bajas son altos y los valores medidos para concentraciones altas son bajos y la región de interés está en el rango de alta concentración.

El cálculo de parámetros admitido por la regresión lineal x/y se utiliza para y.

Se realiza una regresión lineal utilizando los datos transformados.

El resultado se expresa como la pendiente k y la intercepción d.

A partir de este resultado se pueden determinar A = 1/k y B = d/k.

Este cálculo requiere un mínimo de 2 puntos base.

Si fuera necesario, también se puede implementar extrapolación en este proceso. En una evaluación con un grupo experimental, los coeficientes A y B sirven como variables en las transformaciones. Además, también se puede implementar el coeficiente de correlación r. La aproximación siempre se aplica a los valores punto base no escalados.

En el cuadro de diálogo Curva estándar se muestra la información adicional siguiente:

Fórmula con valores de coeficientes A y B

Desviación cuadrática media d

Valor absoluto del coeficiente de correlación r



ADVERTENCIA

ESTA APROXIMACIÓN NO ES CONTINUA Y, POR LO TANTO, UNOS DATOS INCORRECTOS PUEDEN DAR COMO RESULTADO DESVIACIONES ENORMES ENTRE LOS PUNTOS BASE DADOS Y LOS PUNTOS CALCULADOS.

Descripción matemática

Función de aproximación:

$$f:[x_1,x_n] - \{-b\} \to R$$

$$x \mapsto \frac{A \cdot x}{B + x}$$

donde A y B se determinan solucionando el problema de regresión lineal para los puntos base transformados

$$(x_1, \frac{x_1}{y_1}), \dots, (x_n, \frac{x_n}{y_n})$$



Regresión lineal:

$$g:[x_1, x_n] \to R$$

$$x \mapsto k \cdot x + d , \text{ minimizando}$$

$$err(k, d) = \sum_{i=1}^n (g(x_i) - \frac{x_i}{y_i})^2$$

Los parámetros A y B se calculan a partir de k y d según

$$A = \frac{1}{k} \text{ y } B = \frac{d}{k}$$

Esta función f no es continua en –B.

11.4.7 Polinomio

2.º orden: parabólico o cuadrático

3.° orden: cúbico

El usuario puede seleccionar entre el cálculo con un polinomio de orden 2 o el cálculo con un polinomio de orden 3.

Se requiere un mínimo de 3 puntos base para el cálculo con un polinomio de orden 2, mientras que se requiere un mínimo de 4 puntos para el cálculo con un polinomio de orden 3.

Si fuera necesario, en este proceso también se puede implementar extrapolación. En las transformaciones con un polinomio de orden 2 se pueden utilizar los parámetros A = a_0 , B = a_1 y C = a_2 . Para los polinomios de orden tres, se pueden implementar las variables A = a_0 , B = a_1 , C = a_2 y D = a_3 .

En el cuadro de diálogo Curva estándar se muestra la información adicional siguiente:

Fórmula con valores de coeficientes A, B y C para polinomios de orden 2 o fórmula con valores de coeficientes A, B, C y D para polinomios de orden 2

Desviación cuadrática media

Descripción matemática

Función de aproximación:

$$f:[x_1, x_n] \to R$$
$$x \mapsto \sum_{i=0}^{order} a_i \cdot x^i$$
, orden = 2 o 3

(n>orden)

donde aorden,..., ao se determinan minimizando la función de error

$$err(a_{order},...,a_0) = \sum_{i=1}^n (f(x_i) - y_i)^2$$

La solución es única si

$$rank \begin{pmatrix} x_1^{order} & \dots & x_1 & 1 \\ \vdots & \vdots & \vdots & \vdots \\ x_n^{order} & \dots & x_n & 1 \end{pmatrix} = order + 1$$
, que es verdadero si
 $x_i \neq x_j \forall i, j = 1, \dots, n$



11.4.8 Spline cúbica

Los puntos base adyacentes se unirán mediante el cálculo con un polinomio de orden 3. La determinación de los parámetros se lleva a cabo mediante el uso de la condición "not-a-knot".

Este cálculo requiere un mínimo de 3 puntos base.

La extrapolación no es posible. En las transformaciones no hay ningún parámetro que pueda usarse como variable.

Descripción matemática

Características de la función de interpolación:

Polinomio en trozos de orden 3.

Segunda derivación continua en todos los puntos base.

$$\int_{x_1}^{x_n} \left(\frac{d^2 f}{dx^2}\right)^2 dx$$

Condición "not-a-knot": a x2 y xn-1 también la tercera derivación es continua.

11.4.9 Akima

La interpolación Akima crea en algunos casos una curva más suave que la interpolación spline.

Este cálculo requiere un mínimo de 3 puntos base.

La extrapolación no es posible. En las transformaciones no hay ningún parámetro que pueda usarse como variables.

Descripción matemática

Este método de interpolación utiliza polinomios en trozos de orden 3.

El propio polinomio de interpolación en trozos y su primera derivación son continuos.

Método

Las pendientes m_i en todos los puntos base se estiman desde los puntos contiguos por medio de una función especial:

$$q_i = \frac{y_i - y_{i-1}}{y_i - y_{i-1}}$$

 $x_i - x_{i-1}$ son las pendientes del interpolador lineal entre el punto i y el punto i-1, i=2,...,n

$$m_{i} = \frac{q_{i} |q_{i+2} - q_{i+1}| + q_{i+1} |q_{i} - q_{i-1}|}{|q_{i+2} - q_{i+1}| + |q_{i} - q_{i-1}|}, \text{ i=3,...,n-2}$$

Casos especiales:

Para $\mathbf{q}_i = \mathbf{q}_{i+1}$, establecer $\mathbf{p}'(\mathbf{x}_i) = \mathbf{q}_i = \mathbf{q}_{i+1}$.

Para $\mathbf{q}_{i+1} = \mathbf{q}_i$ and $\mathbf{q}_{i+1} \square \mathbf{q}_{i+2}$, resulta $\mathbf{y}_i^* = \mathbf{q}_i$ (analogo para $\mathbf{q}_{i+1} = \mathbf{q}_{i+2}$).

$$m_i = \frac{q_i + q_{i+1}}{2}$$

Para $\mathbf{q}_{i+1} = \mathbf{q}_i$ y $\mathbf{q}_{i+1} = \mathbf{q}_{i+2}$, con

Para los índices 1, 2, n-1, n, este algoritmo no puede estimar las pendientes.



Ahora tenemos las siguientes condiciones para el polinomio de interpolación de tercer orden p_i , i=1, ..., n-1

 $\begin{array}{lll} p_{i}(x_{i}) &= y_{i} \\ p_{i}(x_{i+1}) &= y_{i+1} \\ p_{i}^{\prime}(x_{i}) &= m_{i} \\ p_{i}^{\prime}(x_{i+1}) &= m_{i+1} \end{array}$

que son cuatro condiciones para cada polinomio de interpolación de tercer orden $\ensuremath{p_i}\xspace.$

11.4.10 LogitLog

LogitLog se ha diseñado para curvas estándar sigmoideas. LogitLog siempre producirá una curva con forma de S. La curva alcanzará asintóticamente unos límites superior e inferior, que se pueden establecer independientemente como límites mínimo y máximo. Si el usuario aquí no ha establecido ningún límite, se usarán los valores del punto base mínimo y máximo. La aproximación siempre se aplica a los valores punto base no escalados.

Este cálculo requiere un mínimo de 4 puntos base.

Si fuera necesario, también se puede implementar extrapolación en este proceso. Los parámetros A, B, C y D se pueden emplear como parámetros en las transformaciones.

En el cuadro de diálogo Curva estándar se muestra la información adicional siguiente:

Valores de coeficientes A, B, C y D

Desviación cuadrática media d

Valor absoluto del coeficiente de correlación r (correlación de datos y ajuste)

Restricciones para la aplicación de proceso

LogitLog y la aproximación de Cuatro parámetros solo se pueden emplear al trabajar con curvas estándar que cumplen los siguientes requisitos:

- El proceso sigmoideo aplicado a los datos estándar con respecto a las concentraciones
- Curva generada simétricamente con respecto al valor IC-50
- Para calcular la curva correctamente es necesario tener un valor de concentración de cero y un valor de concentración estándar infinita (valor saturado), de modo que se puedan determinar las asíntotas. Si estos dos valores no están presentes dentro de la curva estándar, se pueden definir manualmente utilizando la opción (Mín, Máx) de Más parámetros de curva estándar. Si no hay valores de Mín. y Máx. definidos, magellan basará el proceso en los valores presentes menor y mayor, respectivamente.



ADVERTENCIA SI NO SE CUMPLEN ESTOS REQUISITOS, UNA APROXIMACIÓN DE ESTA NATURALEZA NO SE PUEDE USAR CORRECTAMENTE.

Descripción matemática

La regresión LogitLog es una función

$$f:[x_1, x_n] \to R$$
$$x \mapsto D + \frac{A - D}{1 + \left(\frac{x}{C}\right)^B}$$

para la descripción de correlación sigmoidea de datos.

El parámetro se puede interpretar como:

$$A = \lim_{x \to 0} f(x)$$
$$D = \lim_{x \to \infty} f(x)$$
$$f(C) = \frac{A+D}{2}$$

A, D se determinan como el mínimo y máximo, respectivamente (o viceversa si la función es decreciente).

Después se resuelve el problema de regresión lineal para los puntos base transformados.

$$X = \log_{10} x$$
$$Y = \ln \left(\frac{\frac{D - y}{D - A}}{1 - \frac{D - y}{D - A}} \right)$$

$$g:[X_i,Y_i]\to R$$

 $x \mapsto k \cdot x + d$, minimizando la función de error

$$err(k,d) = \sum_{i=1}^{n} (g(x_i) - Y_i)^2$$

Los parámetros B, C se determinan a partir de k, d:

$$B = -m \cdot \log_{10}(e)$$
$$C = e^{\frac{b}{B}}$$

Los parámetros A, B, C, D tienen interesantes interpretaciones geométricas.

A = respuesta a una dosis (valor x) de cero (tenga en cuenta que x=0 no se muestra en gráficos logarítmicos)

D = respuesta a una dosis infinitamente alta

C = es el punto de simetría en la curva (IC50); y por debajo de este punto la curva es una imagen reflejada

B = es una medida de la inclinación de la curva en el punto de inflexión; de hecho es (dentro de un posible cambio de signo) la pendiente de la curva escrita en su forma Logit

Tenga en cuenta que la curva siempre es decreciente (si A>D) o creciente (si A<D).



11.4.11 Cuatro parámetros

La aproximación de cuatro parámetros necesita cumplir los mismos requisitos para los puntos base que LogitLog.

La curva resultante se generará utilizando el método de Nelder y Mead Downhill-Simplex. Este método produce una aproximación cada vez más precisa del resultado a lo largo del intervalo. El método de cuatro parámetros requiere considerablemente más tiempo de proceso. Como en el caso de LogitLog, se produce una curva en forma de S.

Este cálculo requiere un mínimo de cuatro puntos base.

Si fuera necesario, también se puede implementar extrapolación en este proceso. Los parámetros A, B, C y D se pueden emplear como parámetros en las transformaciones. La aproximación siempre se aplica a los valores punto base no escalados.

En el cuadro de diálogo Curva estándar se muestra la información adicional siguiente:

Valores de coeficientes A, B, C y D

Desviación cuadrática media d

Valor absoluto del coeficiente de correlación r (correlación de datos y ajuste)

Restricciones para la aplicación de proceso

Para obtener información adicional, consulte la sección 11.4 Tipos de análisis de la curva estándar - LogitLog.

Descripción matemática

Función de aproximación:

$$f:[x_1, x_n] \to R$$
$$x \mapsto D + \frac{A - D}{1 + \left(\frac{x}{C}\right)^B}$$

Método

Primero se calcula una aproximación LogitLog.

Los parámetros A, B, C, D se optimizan con el algoritmo Downhill-Simplex, un algoritmo de minimización que solo necesita evaluaciones de función, ninguna derivación de función.

El algoritmo se detiene si se alcanza (éxito) una precisión de 0,001 o si se supera (fallo) el número máximo de 10.000 iteraciones antes de alcanzar la precisión cada.

Bibliografía

Nelder, J.A., y Mead, R. 1965, Computer Journal, vol. 7, pp 308-313

11.4.12 Cuatro parámetros – Marquardt

La aproximación de cuatro parámetros necesita cumplir los mismos requisitos para los puntos base que LogitLog.

La curva resultante se generará utilizando el método de Levenberg-Marquardt. Este método produce una aproximación cada vez más precisa del resultado a lo largo del intervalo. El método de cuatro parámetros requiere considerablemente más tiempo de proceso. Como con LogitLog, se produce una curva en forma de S.



Este cálculo requiere un mínimo de cuatro puntos base.

Si fuera necesario, también se puede implementar extrapolación en este proceso. Los parámetros A, B, C y D se pueden emplear como parámetros en las transformaciones. La aproximación siempre se aplica a los valores punto base no escalados.

En el cuadro de diálogo Curva estándar se muestra la información adicional siguiente:

- Valores de coeficientes A, B, C y D
- Desviación cuadrática media d
- Valor absoluto del coeficiente de correlación r (correlación de datos y ajuste)

Restricciones para la aplicación de proceso

Para obtener información adicional, consulte la sección 11.4 Tipos de análisis de la curva estándar - LogitLog.

Descripción matemática

Función de aproximación:

$$f:[x_1, x_n] \to R$$
$$x \mapsto D + \frac{A - D}{1 + \left(\frac{x}{C}\right)^B}$$

Método

Primero se calcula una aproximación LogitLog.

Los parámetros A, B, C, D se optimizan con el algoritmo de Levenberg-Marquardt, técnica iterativa que calcula un mínimo local de una función que se expresa como la suma de cuadrados de funciones no lineales.

El algoritmo se detiene si se alcanza (éxito) una precisión de 1E-7 (FLT_EPSILON) o si se supera (fallo) el número máximo de 30.000 iteraciones antes de alcanzar la precisión cada.

11.4.13 Cinco parámetros – Marquardt

La aproximación de cinco parámetros necesita cumplir básicamente los mismos requisitos para los puntos base que LogitLog. No obstante, la curva sigmoidea puede ser no simétrica.

La curva resultante se generará utilizando el método de Levenberg-Marquardt. Este método produce una aproximación cada vez más precisa del resultado a lo largo del intervalo. El método de cinco parámetros requiere considerablemente más tiempo de proceso. Como con LogitLog, se produce una curva en forma de S.

Este cálculo requiere un mínimo de cinco puntos base.

Si fuera necesario, también se puede implementar extrapolación en este proceso. Los parámetros A, B, C, D y E se pueden emplear como parámetros en las transformaciones. La aproximación siempre se aplica a los valores punto base no escalados.

En el cuadro de diálogo Curva estándar se muestra la información adicional siguiente:

- Valores de coeficientes A, B, C, D y E
- Desviación cuadrática media d
- Valor absoluto del coeficiente de correlación r (correlación de datos y ajuste)



Restricciones para la aplicación de proceso

Para obtener información adicional, consulte la sección 11.5 LogitLog.

Descripción matemática

Función de aproximación:

$$f: [x_1, x_n] \to R$$
$$x \mapsto D + \frac{A - D}{(1 + (\frac{x}{C})^B)^E}$$

Método

Los parámetros A, B, C, D, E se optimizan con el algoritmo de Levenberg-Marquardt, técnica iterativa que calcula un mínimo local de una función que se expresa como la suma de cuadrados de funciones no lineales.

El algoritmo se detiene si se alcanza (éxito) una precisión de 1E-7 (FLT_EPSILON) o si se supera (fallo) el número máximo de 30000 iteraciones antes de alcanzar la precisión cada.

11.4.14 Ponderación para ajuste de cuatro / cinco parámetros – ajuste polinomial / Marquardt

La ponderación influye en el algoritmo de optimización estándar.

Dicho algoritmo intenta optimizar la suma ponderada de los mínimos cuadrados.

Un factor de ponderación de 1 significa sin ponderación.

Los factores de ponderación mayores que 1 significan que el punto especificado tiene mayor prioridad para el ajuste resultante.

Los factores de ponderación menores que 1 indican que el punto se tiene menos en cuenta.



ADVERTENCIA

SOLO SE DEBERÁN UTILIZAR LAS PONDERACIONES SI HAY UNA FUERTE EVIDENCIA ESTADÍSTICA Y LOS DATOS RECOPILADOS Y EL ALGORITMO SON APROPIADOS.

Descripción matemática

Desviación cuadrática ponderada media

$$d = \sqrt{\frac{1}{\sum_{i=1}^{n} k_i} \sum_{i=1}^{n} k_i (y_i - f(x_i))^2}$$

Cálculo automático de pesos utilizando varianza:

Los pesos se calculan automáticamente mediante 1/SD², donde SD es la desviación estándar de las réplicas del punto base actual.

Esto significa que los datos con desviación estándar alta tienen menor peso que los datos con desviación estándar baja.



ADVERTENCIA

LA PONDERACIÓN UTILIZANDO VARIANZA ES PELIGROSA CUANDO SE UTILIZAN MUY POCAS RÉPLICAS, YA QUE ENTONCES LA DESVIACIÓN ESTÁNDAR PUEDE TENER UNA PORCIÓN AL AZAR ALTA QUE INFLUYE CONSIDERABLEMENTE EN EL ALGORITMO DE OPTIMIZACIÓN.

Cálculo automático de ponderaciones utilizando ponderaciones relativas:

Los pesos se calculan automáticamente mediante 1/Y², donde Y es el valor medio del punto base actual. Esto significa que el algoritmo minimiza las distancias relativas de los datos a la curva.

11.5 Cálculo de series de dilución

11.5.1 Detección de series de dilución

Es preciso cumplir los siguientes criterios para forzar a **magellan** a detectar series de dilución en el diseño:

- Muestra(s) con un mínimo de cuatro réplicas
- Utilización de al menos cuatro factores de dilución distintos para las réplicas simples de una muestra

magellan comprueba el diseño completo y selecciona todas las series de dilución encontradas.

11.5.2 Cálculo de parámetros de curvas

Los parámetros de las curvas se calculan para cada serie de dilución encontrada utilizando el algoritmo **Cuatro parámetros – Marquardt**. En caso de fallo, se vuelve a procesar el cálculo utilizando el algoritmo **Cuatro parámetros**. Consulte en la sección 11.4 Tipos de análisis de la curva estándar información detallada acerca de los algoritmos mencionados anteriormente.

11.5.3 Cálculo de valores de IC

Primero se seleccionan la intensidad máxima $(I_{\text{máx}})$ y la intensidad mínima $(I_{\text{mín}})$ de cada serie de dilución.

Cálculo de intensidad IC

Si el valor de 0% se ajusta a intensidad "Intensidad 0"

$$I_{IC} = \frac{ICx}{100.0} * I_{\max}$$

Si el valor de 0% se ajusta a "Intensidad mín. de las series de dilución"

$$I_{IC} = I_{\min} + \frac{(I_{\max} - I_{\min}) * ICx}{100.0}$$

Cálculo de dilución a ICx

La dilución resultante se calcula utilizando los parámetros de las curvas calculados en el paso 12.6.2 y la intensidad IC calculada.

dilution =
$$C * (\frac{A - I_{IC}}{I_{IC} - D})^{\frac{1}{B}}$$

12. Ejemplo de aplicación

12.1 Ejemplo paso a paso: ELISA cuantitativo

Este capítulo contiene un ejemplo paso a paso (prueba cuantitativa) de cómo crear un método en **magellan**. Siga las instrucciones para aprender a definir evaluaciones a partir de la descripción de un kit de prueba en **magellan**.

El método se creó con un Sunrise en modo demo, pero también se pueden definir los métodos con otro instrumento Tecan conectado, siempre que sea capaz de medir la absorbencia.

12.1.1 Descripción del kit de prueba

La descripción del fabricante que acompaña al kit de prueba de un ELISA IgM (detección de anticuerpos) cuantitativo contiene las siguientes instrucciones: Diseño de placa:

1	2	3	4	5	6	7	8	9	10	11	12
BLK	C3	S1									
NC	C4	S2									
NC	C4	S2									
C1	C5	S3									
C1	C5	S3									
C2	C6										
C2	C6										
C3	S1										
	1 BLK NC C1 C1 C2 C2 C2 C3	1 2 BLK C3 NC C4 NC C4 C1 C5 C1 C5 C2 C6 C3 S1	1 2 3 BLK C3 S1 NC C4 S2 NC C4 S2 C1 C5 S3 C2 C6 C3 S1	1 2 3 4 BLK C3 S1 NC C4 S2 NC C4 S2 C1 C5 S3 C2 C6 C3 S1	1 2 3 4 5 BLK C3 S1 NC C4 S2 NC C4 S2 C1 C5 S3 C2 C6 C3 S1	1 2 3 4 5 6 BLK C3 S1 NC C4 S2 NC C4 S2 NC C4 S2 C1 C5 S3 C2 C6 C3 S1	1 2 3 4 5 6 7 BLK C3 S1 NC C4 S2 NC C4 S2 NC C4 S2 <td>1 2 3 4 5 6 7 8 BLK C3 S1 <</td> <td>1 2 3 4 5 6 7 8 9 BLK C3 S1 NC C4 S2 NC C4 S2 NC C4 S2 NC C4 S2 </td> <td>1 2 3 4 5 6 7 8 9 10 BLK C3 S1</td> <td>1 2 3 4 5 6 7 8 9 10 11 BLK C3 S1 <</td>	1 2 3 4 5 6 7 8 BLK C3 S1 <	1 2 3 4 5 6 7 8 9 BLK C3 S1 NC C4 S2 NC C4 S2 NC C4 S2 NC C4 S2 	1 2 3 4 5 6 7 8 9 10 BLK C3 S1	1 2 3 4 5 6 7 8 9 10 11 BLK C3 S1 <

BLK = Blanco, NC = Control negativo, C1 - C6 = Calibradores (Estándares),

S1 – S... = Muestras

Medición y evaluación

Leer placa a una longitud de onda de 492 nm, referencia a 620 nm.

Placa/lector del blanco en pocillo A1.

Concentraciones de los calibradores (Estándares):

Calibrador 1	5 UA/mL
Calibrador 2	10 UA/mL
Calibrador 3	20 UA/mL
Calibrador 4	40 UA/mL
Calibrador 5	80 UA/mL
Calibrador 6	160 UA/mL

Después de la corrección del blanco se trazan las densidades ópticas (DO492– DO620) como función de la concentración. La línea de regresión que atraviesa estos puntos es la curva estándar. Interpretación de los resultados de la prueba:

IgM < 18 UA/mL	Negativo
18 UA/mL <= IgM < 22 UA/mL	Intermedio
IgM >= 22 UA/mL	Positivo

La concentración de IgM calculada de ambos controles negativos debe ser inferior a 8 UA/mL.

Procesamiento de datos

Después de la medición, el archivo de datos (zona de trabajo) se guarda automáticamente y se crea un informe que contiene los parámetros de medición, diseño de placa, valores de blancos, curva estándar, concentraciones de IgM, definición de corte, resultados cualitativos de las muestras y validaciones.

Además, la distribución y los resultados cualitativos se guardarán en un archivo ASCII.

12.1.2 Crear un método

En el cuadro de diálogo Lista de Asistentes seleccione Crear o editar método y haga clic en Aceptar. En la página Bienvenida del Asistente Crear o editar método haga clic en Siguiente y se mostrará el cuadro de diálogo Seleccionar un archivo. Elija Crear nuevo.

Parámetros de medición

Haga clic en Siguiente para ver el cuadro de diálogo Parámetros de medición.

Crear o editar métod	lo	×
	Tipo de m	edición Medición de punto final Medición cinética Medición cinética de pocillo Medición de múltiples etiquetas Parámetros de medición
Ayuda	<<< Atrás	DEFINIR EVALUACIÓN
		© 2016 Tecan

Seleccione **Medición de punto final** y haga clic en el botón **Parámetros de medición**. Se muestra el siguiente cuadro de diálogo:



arámetro	de medición (SUNF	RISE Simulatio	n)			×
General	Longitudes de onda	Parámetros d	e me <mark>dic</mark> ión	Temperatura	Agitación	
) Longitud de onda fija					
С) Exploración de absor	bencia	Establ	lecer rango de onda comp	longitudes de leto	
Longi	tud de onda					
Lor Lor	ngitud de onda <mark>d</mark> e med ngitud de onda de refe	lición: renc <mark>i</mark> a:	492 620	nm nm		
					OK Canad	Hab

En la pestaña General, seleccione Absorbencia.

En la pestaña **Longitudes de onda**, seleccione 492 nm como Longitud de onda de la medición y 620 nm como Longitud de onda de referencia.

Haga clic en **Aceptar** para regresar al cuadro de diálogo **Parámetro de medición**.

Haga clic en **Definir evaluación** y se mostrará la ventana **Diseño de las placas**.





Configurar diseño

Defina el diseño de placa en el cuadro de diálogo **Asignación de pocillos** del lado derecho de la pantalla.

En el cuadro de grupo Identificadores, seleccione BL (Blanco).

En el cuadro de grupo Experimental permanece el número 1.

En el cuadro de grupo Réplicas, seleccione Todas.



Haga clic en el pocillo A1, que quedará marcado con un borde rojo.

Haga clic en **Rellenar selección** y el pocillo quedará marcado con el tipo de identificador elegido.



Aviso También se puede llenar por separado un pocillo si se hace doble clic sobre el mismo.

A continuación, elija los siguientes ajustes en el cuadro de diálogo **Asignación** de pocillos:

En el cuadro de grupo Identificadores, seleccione NC (Control Negativo).

En el cuadro de grupo Experimental permanece el número 1.

Se selecciona Todos en el cuadro de grupo Réplicas.

Empezando en el pocillo **B1** haga clic y arrastre el ratón hasta **C1**. Los pocillos **B1** a **C1** quedarán marcados con un borde rojo.

Haga clic en **Rellenar selección** y los pocillos quedarán marcados con el tipo de identificador elegido.

Luego se deben asignar los calibradores (estándares) a los pocillos **D1** a **G2**. Seleccione los siguientes ajustes del cuadro de diálogo **Asignación de pocillos**:

En el cuadro de grupo Identificadores, seleccione ST (Estándar).

En el cuadro de grupo Experimental permanece el número 1.



En el cuadro de grupo Réplicas, seleccione entre Número fijo y Todas:

Número fijo

Aplicable solamente a estándares y muestras que admiten el uso de IDs. Si está activo el botón **Número fijo**, se puede escribir un número en el campo de texto correspondiente. Ese número define la cantidad de réplicas previstas para el método. En los pocillos seleccionados, se creará el número indicado de réplicas para cada ID. Por lo tanto, el número de pocillos seleccionados deberá ser un múltiplo del número de réplicas que se haya escrito.

Todas

Todos los pocillos seleccionados estarán definidos como réplicas. Si se elige un número de ID existente para muestras y estándares, los pocillos seleccionados se añadirán como réplicas a las réplicas existentes. Con el resto de tipos de identificadores, los pocillos seleccionados se añadirán como réplicas a las réplicas existentes.

Los dos botones con flechas definen la dirección de secuencia (horizontal o vertical) de réplicas y números de ID.

En el presente ejemplo, seleccione Número fijo y 2.

En la casilla **Número-ID** y en el cuadro de grupo **Réplicas** seleccione las **flechas verticales**.

Después seleccione los pocillos D1 a G2 y haga clic en Rellenar selección.



Aviso

Seleccione los pocillos como sigue: empezando en el pocillo D1, haga clic y arrastre el ratón sobre los pocillos requeridos hasta H1. Después mantenga presionada la tecla Control (Ctrl) y arrastre el ratón sobre los pocillos requeridos de A2 a G2.

El Diseño de placa aparecerá así:

Create/Edit a Method														
Undo Redo Select all unu	used	10%	100%) 🗆 Zoom										
(()	0													
Method layout Image: Constraint of the second	eate/edit n		1	2	3	4	5	6	7	8	9	10	Well Assignm. Identifiers	Exp. group:
Transformed data 🛞	ethod	A	BL1 1/1	ST1_3 2/2									SM BL ST	ID-Num.
Concentrations 🔗		B	NC1 1/2	ST1_4 1/2									PC NC	
Conc. transformed d (*) Add new conc. transformati		C	NC1 2/2	ST1_4 2/2									LPC HPC	Replicates →
Evaluate data 🛞		D	ST1_1 1/2	ST1_5 1/2									BF RF	Fix number:
QC Validation Data handling		E	ST1_1 2/2	ST1_5 2/2									Fill selection	Delete
Data export Printed report Automated data handling		F	ST1_2 1/2	ST1_6 1/2										
Miscellaneous 📀 User prompts		G	ST1_2 2/2	ST1_6 2/2										CAN Austri
Vumber format Method notes		Н	ST1_3 1/2											TEC
Help Cancel <<< Back													FINISH	
													@ 2004 Tecap	

2023-07



Para seleccionar todos los pocillos vacíos de la placa, en la barra de herramientas haga clic en **Seleccionar no usados**. A continuación mantenga presionada la tecla Control (Ctrl) y haga clic en el pocillo **H12**, de modo que quede en blanco y sin marcar.

En el cuadro de diálogo **Asignación de pocillos** seleccione **SM (Muestra)** bajo **Identificadores**.

En el cuadro de grupo Experimental permanece el número 1.

En el cuadro de grupo Réplicas seleccione entre Número fijo y 2.

En la casilla **Número-ID** deje 1 y en el cuadro de grupo **Réplicas** seleccione las **flechas verticales**. A continuación haga clic en **Rellenar selección**. El procedimiento de Definición de diseño queda entonces completado.

Transformaciones

Para definir la reducción del blanco, en la barra de control a la izquierda de la ventana seleccione la siguiente opción, **Añadir nueva transformación...** en el elemento **Datos transformados**.



Se muestra un cuadro de diálogo para preguntarle si desea definir una reducción del blanco. Haga clic en **Sí**.

Seleccione **Datos de diferencia** en el recuadro **Entrada**. De forma predeterminada la transformación se denomina Transformation1 (véase barra de control). Si ya ha confirmado antes la definición de una reducción del blanco, el software la llamará **Reducción del blanco** automáticamente.

En el cuadro **Fórmula** se mostrará automáticamente **x-BL1** para esta reducción del blanco, donde x se refiere al valor actual de los datos de entrada en un pocillo y BL1 es el valor medio de los pocillos en blanco del grupo experimental 1.

Ahora haga clic en el signo que hay en la esquina superior izquierda de la microplaca para seleccionar la placa completa y haga clic en la marca de verificación verde junto a la ventana de fórmulas para asignar las transformaciones a los pocillos. Consulte detalles y explicaciones adicionales acerca de la definición y la asignación de transformaciones en la sección 4.3.5 Datos transformados: Añadir nueva transformación. Se muestra la siguiente ventana:



nottado 2 Privado Datos de dife	rencia		¢									~ [Da	tos deponibles (corp	Ce ritse de detos * Tu
pikos anc., DL., Ref demades	0	1	2	3	4	5	6	7	8	9	10	11	12	
States 2	A	BL1 1/1 x-BL1	ST1_3 2/2 x-BL1	SM1_1 2/2 x-BL1	SM1_5 2/2 x-BL1	SM1_9 2/2 x-BL1	SM1_13 2/2 x-BL1	SM1_17 2/2 x-BL1	SM1_21 2/2 x-BL1	SM1_25 2/2 x-BL1	SM1_29 2/2 x-BL1	SM1_33 2/2 x-BL1	SM1_37 2/2 x-BL1	
ter de de la contra		NC1 1/2 x-8L1	ST1_4 1/2 x-BL1	SM1_2 1/2 x-BL1	SM1_6 1/2 x-BL1	SM1_10 1/2 x-BL1	SM1_14 1/2 x-BL1	SM1_18 1/2 x-BL1	SM1_22 1/2 x-BL1	SM1_26 1/2 x-BL1	SM1_30 1/2 x-BL1	SM1_34 1/2 x-BL1	SM1_38 1/2 x-BL1	
is autoratinal	C	NC1 2/2 x-BL1	ST1_4 2/2 x-BL1	SM1_2 2/2 x-BL1	SM1_6 2/2 x-BL1	SM1_10 2/2 x-BL1	SM1_14 2/2 x-BL1	SM1_18 2/2 x-BL1	SM1_22 2/2 x-BL1	SM1_26 2/2 x-BL1	SM1_30 2/2 x-BL1	SM1_34 2/2 x-BL1	SM1_38 2/2 x-BL1	
*		ST1_1 1/2 x-BL1	ST1_5 1/2 x-BL1	SM1_3 1/2 x-BL1	SM1_7 1/2 x-BL1	SM1_11 1/2 x-BL1	SM1_15 1/2 x-BL1	SM1_19 1/2 x-BL1	SM1_23 1/2 x-BL1	SM1_27 1/2 x-BL1	SM1_31 1/2 x-BL1	SM1_35 1/2 x-BL1	SM1_39 1/2 x-BL1	
		ST1_1 2/2 x-BL1	ST1_5 2/2 x-BL1	SM1_3 2/2 x-BL1	SM1_7 2/2 x-BL1	SM1_11 2/2 x-BL1	SM1_15 2/2 x-BL1	SM1_19 2/2 x-BL1	SM1_23 2/2 x-BL1	SM1_27 2/2 x-BL1	SM1_31 2/2 x-BL1	SM1_35 2/2 x-BL1	SM1_39 2/2 x-BL1	
	F	ST1_2 1/2 x-BL1	ST1_6 1/2 x-BL1	SM1_4 1/2 x-BL1	SM1_8 1/2 x-BL1	SM1_12 1/2 x-BL1	SM1_16 1/2 x-BL1	SM1_20 1/2 x-BL1	SM1_24 1/2 x-BL1	SM1_28 1/2 x-BL1	SM1_32 1/2 x-BL1	SM1_36 1/2 x-BL1	SM1_40 1/2 x+BL1	
		ST1_2 2/2 x-8L1	ST1_6 2/2 x-BL1	SM1_4 2/2 x-BL1	SM1_8 2/2 x-BL1	SM1_12 2/2 x-BL1	SM1_16 2/2 x-BL1	SM1_20 2/2 x-BL1	SM1_24 2/2 x-BL1	SM1_28 2/2 x-BL1	SM1_32 2/2 x-BL1	SM1_36 2/2 x-BL1	SM1_40 2/2 x-BL1	AN Austria
	H	ST1_3 1/2 x-BL1	SM1_1 1/2 x-BL1	SM1_5 1/2 x-BL1	SM1_9 1/2 x-BL1	SM1_13 1/2 x-BL1	SM1_17 1/2 x-BL1	SM1_21 1/2 x-BL1	SM1_25 1/2 x-BL1	SM1_29 1/2 x-BL1	SM1_33 1/2 x-BL1	SM1_37 1/2 x-BL1		TEC

En cada pocillo se mostrará la siguiente información (pocillo de ejemplo A5):

SM1_9	Muestra, grupo experimental número 1, ID de muestra número 4.
2/2	El número de réplica es 2, el número total de réplicas es 2.
x-BL1 o 1000	Transformación asignada x-BL1 (cuando se ha seleccionado Transformación) o valor del Factor de dilución de 1 (cuando se han seleccionado Valores de Conc., Dil., Ref.).

Definición de valor de concentración / dilución / referencia

Para definir los valores respectivos como se describe en el kit de prueba, en el elemento **Diseño de método** de la barra de control, seleccione **Valores de Conc., Dil., Ref.**.

Calibrador 1	5 UA/mL
Calibrador 2	10 UA/mL
Calibrador 3	20 UA/mL
Calibrador 4	40 UA/mL
Calibrador 5	80 UA/mL
Calibrador 6	160 UA/mL

Asegúrese de que esté seleccionado ST en la lista Seleccionar identificador.

En la lista **Identificador** se mostrará una lista de los estándares del Grupo exp. 1. En el cuadro correspondiente a **Concentración** de **ST1_1**, escriba el número **5** y en el cuadro **Unidad**, escriba UA/mL. En el cuadro correspondiente a **Concentración** de **ST1_2** escriba el número **10**. La unidad solamente se debe definir una vez y será válida para todos los estándares. Del mismo modo, escriba los valores para ST1_3 a ST1_6.



Image: Second	0	1	2	3	4	5	6	7	8	9	10	11	12
All 40 5 00 reacts banformasi mitraciones	A	BL1 1/1 1	ST1_3 2/2 20	SM1_1 2/2 1	SM1_5 2/2 1	SM1_9 2/2 1	SM1_13 2/2 1	SM1_17 2/2 1	SM1_21 2/2 1	SM1_25 2/2 1	SM1_29 2/2 1	SM1_33 2/2 1	SM1_37 2/2 1
n de conc. Instru-, ;; r nueve bandormado unr datum		NC1 1/2 1		SM1_2 1/2 1	SM1_6 1/2 1	SM1_10 1/2 1	SM1_14 1/2 1	SM1_18 1/2 1	SM1_22 1/2 1	SM1_26 1/2 1	SM1_30 1/2 1	SM1_34 1/2 1	SM1_38 1/2 1
ada QC maarmanta da da <u>con</u> toda de dato na ingenes		NC1 2/2 1	ST1_4 2/2 40	SM1_2 2/2 1	SM1_6 2/2 1	SM1_10 2/2 1	SM1_14 2/2 1	SM1_18 2/2 1	SM1_22 2/2 1	SM1_26 2/2 1	SM1_30 2/2 1	SM1_34 2/2 1	SM1_38 2/2 1
samento automotizad	D	ST1_1 1/2 5	ST1_5 1/2 80	SM1_3 1/2 1	SM1_7 1/2 1	SM1_11 1/2 1	SM1_15 1/2 1	SM1_19 1/2 1	SM1_23 1/2 1	SM1_27 1/2 1	SM1_31 1/2 1	SM1_35 1/2 1	SM1_39 1/2 1
Cristed: Unite:			ST1_5 2/2 80	SM1_3 2/2 1	SM1_7 2/2 1	SM1_11 2/2 1	SM1_15 2/2 1	SM1_19 2/2 1	SM1_23 2/2 1	SM1_27 2/2 1	SM1_31 2/2 1	SM1_35 2/2 1	SM1_39 2/2 1
- s(p+1) = s(p) = 0 - s(p+1) = s(p) = 0 - s(p+1) = s(p) = 0 - s(p+1) =			ST1_6 1/2 160	SM1_4 1/2 1	SM1_8 1/2 1	SM1_12 1/2 1	SM1_16 1/2 1	SM1_20 1/2 1	SM1_24 1/2 1	SM1_28 1/2 1	SM1_32 1/2 1	SM1_36 1/2 1	SM1_40 1/2 1
			ST1_6 2/2 160	SM1_4 2/2 1	SM1_8 2/2 1	SM1_12 2/2 1	SM1_16 2/2 1	SM1_20 2/2 1	SM1_24 2/2 1	SM1_28 2/2 1	SM1_32 2/2 1	SM1_36 2/2 1	SM1_40 2/2 1
	H		SM1_1 1/2 1	SM1_5 1/2 1	SM1_9 1/2 1	SM1_13 1/2 1	SM1_17 1/2 1	SM1_21 1/2 1	SM1_25 1/2 1	SM1_29 1/2 1	SM1_33 1/2 1	SM1_37 1/2 1	

Aparecerá la pantalla que muestra el diseño de placa y la concentración:

Curva estándar

Para definir la curva estándar apropiada, en la barra de control haga clic en **Curva estándar** del elemento **Concentraciones**.

La descripción del kit de prueba indica:

Después de la corrección del blanco, las densidades ópticas (DO492 – DO620) se representan gráficamente frente a la concentración. La línea de regresión que atraviesa estos puntos es la curva estándar.

En la pestaña **Datos**, seleccione **Reducción del blanco** como datos de entrada.

Jurva	estantidar				
atos	Tipo de análisis	Intercepciones	Eje	Gráfico	
	Entrada:	Reducción del bla	anco	<u> </u>	
	_				
	Estándares e	n diseño			
	O Estándares e	n archivo ext:		Selectionar	
	O Estándares e	n grupo experime	ntal:		
	O Sin gráfico de	curva estándar			
	Concentracion	nes adicionales			



Datos	Tipo de análisis Intercepciones Eje Gráfi	0
	Tipo de análisis Punto a punto Regresión lineal Regresión no lineal Estría cúbica Akima	Escala de datos: Lin(x)Lin(y) ~
	Polinómica Cuatro parámetros Marquardt cuatro parámetros Cinco parámetros LogitLog	€ Más
	☐ Induir (0,0) ☐ Factor de extrapolación: 1	

En la pestaña Tipo de análisis, seleccione Regresión lineal.

En la pestaña **Eje**, defina la etiqueta y la escala de los ejes como se muestra a continuación:

Eje X		
Etiqueta: Concentración [U	/mL]	
Color:	Escala log	
Autoseleccionar rango		
◯ Intervalo	mín.: máx.:	
Retícula	Color: Estilo:	
Eje Y		
Etiqueta: Reducción del bla	 C0	
Color:	Escala log	
Autoseleccionar rango		
◯ Intervalo	mín.: máx.:	
Retícula	Color: Estilo:	



En la pestaña **Gráfico**, defina el título del gráfico, las curvas, el tipo de letra y la visualización del gráfico.

atos	Tipo de análisis Intercepciones	Eje Grafico		
	Título			
	Etiqueta: IgM ELISA			
	Color:			
	Curvas			
	Etiq	ueta: Grp. 1		
	Color: Símbolo: Al Tamaño: Medio V	Esconder curva		
	Fuente	Visualizar		
	Pequeño	✓ Leyenda	✓ Intercepciones	
	O Medio O Grande	Puntos base	Barras de errores	

Definir cortes

Para definir los límites para la evaluación cualitativa, en la barra de control seleccione **Definición del corte** en el elemento **Evaluar datos**.

La descripción del kit de prueba contiene las siguientes instrucciones:

Interpretación de los resultados de la prueba:

lgM < 18 UA/mL	Negativo
18 UA/mL <= IgM < 22 UA/mL	Intermedio
IgM >= 22 UA/mL	Positivo

Aplique el siguiente procedimiento para definir los cortes adecuados:

En el cuadro Datos de entrada, seleccione Conc. media (UA/mL).

La tabla **Cortes** representa una escala que indica el extremo alto y bajo para los **Límites** y las **Etiquetas**. En **Límites**, escriba 22 como primer límite (máximo) y 18 como segundo límite (mínimo).

En **Etiquetas**, escriba la interpretación de la prueba (**Positivo**, **Intermedio** y **Negativo**) en los cuadros individuales. Para asignar un color, utilice la paleta de colores desplegable:

Positivo - Rojo

Intermedio – Azul

Negativo – Verde



La pantalla contiene lo siguiente:

itrada:	Conc. media. (UA/mL)		~		
ortes						
Colorea	Etiquetas	Lími	tec			
COIOTES	Positive	≜ ∰	les			
	Intermedia	- 22				
	Negative	18				
		┾┙╉╸┝				
		+-+-				
		┲┹				_
						_
						_
	•	- Tow				
		17				
ntrada de	fórmula				ſ	Prueba competitiva
Variable		Operadore	es Funciones			
BL1	~	+	✓ and	~		

Haga clic en **Seleccionar resultados de corte** para elegir los tipos de identificadores que indicarán los resultados de corte.

Definir validaciones del control de calidad (QC)

En la barra de control, haga clic en **Validación QC** en el elemento **Evaluar datos**. Se deben definir criterios de validación para la prueba de modo que se garantice la validez de los resultados de la prueba.

En el ejemplo dado se deben cumplir los siguientes requisitos:

La concentración de IgM calculada de ambos controles negativos debe ser inferior a 8 UA/mL.

En el cuadro Entrada, seleccione Conc. simple (UA/mL).

En la primera fila escriba NC1_1<8 o escriba una fórmula usando las variables, operadores y funciones disponibles.



Aviso NC1_1 significa Control negativo del grupo experimental 1, réplica 1.

En la segunda fila, escriba NC1_2<8.



12. Ejemplo de aplicación

Entonces se mostrará el cuadro de diálogo Validación QC como sigue:

	Entrada:	Cond	: , simple	. (UA/r	nL)			~		Grupo	<mark>d</mark> e valida	ación: [1
	Condicio	es de	validad	ión		 				 			
	NC1_1<8												
Ś	NC1_2<8												
N.													
ŝ													
5													
ŝ										 			
2						 	 		 				
	I												

Organizar el informe impreso

En la barra de control, haga clic en **Informe impreso** del elemento **Procesamiento de datos**. Se muestra la pantalla siguiente:

Imprimir como O Lista	Puede utilizar arrastrar y colocar para crear un informe. Seleccione un elemento en la lista de datos disponible y arrástrelo a la lista de datos seleccionada. Colóquelo en la posición apropiada dentro del informe.
Datos disponibles:	Datos seleccionados:
Datos del instrumento Datos del instrumento Datos reducidos Datos de diferencia Datos de diferencia - Media Datos de diferencia - Desviac Datos de diferencia - Coeficie	Insertar Anexar Anexar Anexar Definition Definition Definition Definition Definition Definition Definition Definition Definition Definition Definition Definition Definition Definition Definition Definition Definition Definition Definition Definition Definition Definition Definition Definition Definition Definition Definition Definition Definition Definition Definition Definition Definition Definition Definition Definition Definition Definition Definition Definition Definition Definition Definition Definition Definition Definition Definition Definition Definition Definition Definition Definition Definition Definition Definition Definition Definition Definition Definition Definition Definition Definition Definition Definition Definition Definition Definition Definition Definition Definition Definition Definition Definition Definition Definition Definition Definition Definition Definition Definition Definition Definition Definition Definition Definition Definition Definition Definition Definition Definition Definition Definition Definition Definition Definition Definition Definition Definition Definition Definition Definition Definition Definition Definition Definition Definition Definition Definition Definition Definition Definition Definition Definition Definition D
	Eliminar Eliminar todos
R Notas E Protocolo de errores E Registro de modificaciones E Firma	

En la pestaña **Selección de datos**, el cuadro **Datos disponibles** contiene todos los datos de informe disponibles. Con los botones **Insertar y Anexar**, transfiera datos al cuadro **Datos seleccionados**. Para transferir datos, también puede arrastrar y colocar.

En la casilla **Imprimir como**, seleccione si imprimirá los datos como una matriz o como una lista con una orientación especial.

En el ejemplo dado se debe crear un informe que contenga los parámetros de medición, el diseño de placa, valores de blanco, curva estándar, concentraciones IgM, definición de corte, resultados cualitativos de las muestras y validaciones.

Antes de crear el informe, deberá quitar **Matriz-Datos de diferencia** que aparece de modo predeterminado en la casilla **Datos seleccionados**. Solamente **Parámetros de medición** permanece en el cuadro **Datos seleccionados**.



Seleccione **Diseño de método/Diseño** en el cuadro **Datos disponibles** y adjúntelo como una matriz al informe. Para ello, haga clic en **Anexar**. Después, inserte **Reducción del blanco**, **Conc. media (UA/mL)** y **Resultados de corte** en la matriz. Para ello, seleccione los elementos correspondientes y haga clic en **Insertar**.

Añada Gráfico: Curva estándar, Definición de corte y Criterios de validación QC a los datos seleccionados. La parte de configurar datos del procedimiento de definición del informe está lista; el cuadro de diálogo Informe impreso tiene este aspecto:



En las pestañas **Encabezado** y **Pie de página** defina el diseño del encabezado y del pie de página del informe (consulte los detalles en la sección 4.3.13 Procesamiento de datos: Informe impreso).

Exportar datos

En la barra de control, seleccione **Exportación de datos** en el elemento **Procesamiento de datos**. En este ejemplo, el diseño y los resultados de corte se deberán guardar como un archivo ASCII. Seleccione **Diseño** y **Resultados de corte** en la ventana **Datos disponibles**; haga clic en la flecha \rightarrow para insertarlos en la ventana **Datos seleccionados**. La pantalla muestra la siguiente información:



1 N. 2	-		
Datos reducidos Datos de diferencia Datos de diferencia - Media Datos de diferencia - Desviación e Datos de diferencia - Coeficiente e Datos transformados Datos transformados Concentraciones Resultados de corte ', ID de muestra Diseño de método Diseño de métodos Resultaciones de los pocillos Nombres para métodos de banda D Concentraciones originales Factores de dilución Diseño	s Je	-> <-	abajo
< >			



Aviso Los datos exportados siempre deberán contener la lista IDs de muestras o Diseño.

Procesamiento automatizado de datos

En la barra de control, seleccione **Procesamiento automatizado de datos** en el elemento **Procesamiento de datos**.



Seleccione **Exportar a archivo ASCII** y Ver resultados después de la **medición**. En magellan **Tracker, Guardar zona de trabajo** está seleccionado de modo predeterminado y no se puede modificar.



Guardar el método

Haga clic en **FINALIZAR** para abrir la ventana **Guardar como**. Escriba el nombre de archivo del método y, si fuera necesario, complete los demás campos.

Create/Edit a Method				×				
Save in: mth Filename: Application Example_en.mt File remarks: Audit trail comment: Organize favorites Audit t Helo	Name Application Example_ch Application Example_de Application Example_es Application Example_is Application Example_ja Application Example_po Application Example_po Application Example_ru Cx Cx_SUNRISE Multilabel_SUNRISE Spectrum_SUNRISE Spectrum_SUNRISE	Remarks	Status Unsigned Unsigned Unsigned Unsigned Unsigned Unsigned Unsigned Unsigned Unsigned Unsigned Unsigned Unsigned					
Cancel <<< Back	Run	this method now	SAVE&FINISH					
Campo de texto Nombre de archivo	Se debe escribir un nor automáticamente un no predeterminado, pero s	mbre de arc ombre de ar se puede ca	chivo. Se su rchivo ambiar.	giere				
Campo de texto Comentarios del archivo	Las observaciones aqu mostrarán con el nomb	uí anotadas pre del archi	se guardará vo.	àn y				
Campo de texto Comentarios al registro de modificaciones	Los comentarios aquí anotados serán almacenados en el registro de modificaciones. <i>Esta opción solamente está disponible con</i> magellan Tracker.							
Botón Registro de modificaciones	Se muestra el cuadro o modificaciones. Esta con magellan Tracker	de diálogo F opción sola	Registro de mente está	disponible				
Botón Organizar favoritos	Se muestra el cuadro de diálogo Organizar favoritos . (Consulte la sección 6.4 Activar favorito – Organizar favoritos).							
Botón Firmas	Se muestra el cuadro o solamente está dispon	de diálogo F <i>ible con</i> ma	Firma. Esta gellan Trac	opción : ker .				
Contraseña de método	Introduzca una contras proteger el método que sección 4.4.1 Protecció contraseña).	eña para es e va a guarc ón de métoc	se método s dar (consulte dos mediante	i desea e la e				
Casilla de verificación Ejecutar este método ahora	El método se ejecutará finalice el asistente.	inmediatar	mente despu	ués de que				



Registro de modificaciones

Haga clic en **Registro de modificaciones** para abrir el cuadro de diálogo **Registro de modificaciones**:

	Usuario	Fecha	Acción	Comentario	
1	Administrator (Admin)	2018-03-06 08:52:42	Creado		
2	Administrator (Admin)	2018-03-06 09:59:10	Modificado		
3	Administrator (Admin)	2018-03-07 09:07:10	Modificado		

El registro de modificaciones muestra una lista de todas las modificaciones del método efectuadas. Cada anotación consta de los datos de usuario (nombre y nombre completo), fecha y hora de la modificación, si el archivo ha sido creado o modificado, así como comentarios al registro de modificaciones.

Haga clic en **Presentación preliminar...** para acceder a una vista previa del archivo. Si desea comparar un método con su versión anterior deberá primeramente imprimir esta, ya que no es posible abrir dos vistas preliminares de impresión simultáneamente.

12.1.3 Ejecutar el método

Si está seleccionado **Ejecutar este método ahora** en el cuadro de diálogo **Guardar como** del **Asistente Crear o editar método**, se mostrará el cuadro de diálogo **Asistente Iniciar medición/Iniciar medición** después de hacer clic en **Guardar**:

Medicion		Parámetros de medición			
Zona de trabajo: 07032018-008.wsp		Ciclo cin. arb Modo de medición: Absorbencia Longitud de onda de medición: 492 nm			
vietodo: ista da IDa da musatra:	Application Example_es.mt	Modificar diseño	Longitud de onda de referencia: 620 nm		
usta de ilos de muestra:		Insertar	Modo lectura: normal Unidad: OD		
nstrumento	/				
Utilizar apilador		Actual: 32.1 °C			
Introducir placa	Control de temperatura	Objetivo: 37.0 °C			
Movimientos	Control inyector	Optimiza la posición Z			
Tenga en cuenta lo sigui	ente:				
	/				
Ayuda					


Para empezar la medición, haga clic en **Iniciar**. Se creará automáticamente una zona de trabajo que contiene toda la información previamente introducida y que recogerá todos los valores de medición. Durante la ejecución de la medición se mostrará un cuadro de diálogo de estado de la medición que indica el progreso de la misma.

Tras completarse la medición, aparecerá el cuadro de diálogo **Resultados** donde se muestran los resultados y los cálculos.

12.1.4 Evaluar el resultado

Evaluar resultados le permite ver y evaluar datos sin procesar. Se pueden visualizar los parámetros de evaluación, así como volver a evaluar los datos.

Esta sección le guía a través del **asistente Evaluar resultados** usando dos archivos de zonas de trabajo de ejemplo creados ejecutando el método ELISA cuantitativo.



Aviso

En la Lista de métodos de magellan Standard se muestran automáticamente archivos de ejemplo. Para magellan Tracker, esos archivos están disponibles en el directorio de datos predeterminado y se deben convertir.

En el cuadro de diálogo Lista de asistentes, haga clic en Evaluar resultados.

Haga clic en Siguiente en la página Bienvenida del Asistente Evaluar resultados y se mostrará el cuadro de diálogo Seleccionar un archivo.

Seleccione la zona de trabajo **Quantitative Elisa example_Sunrise.wsp** en la lista de archivos y haga clic en **Haga su selección**. Se ejecutarán los cálculos y se mostrará la siguiente ventana de diseño de placa:



En cada pocillo individual se mostrará el valor calculado. En función del elemento seleccionado en la barra de control, la ventana de diseño de placa cambiará como corresponda. Los parámetros y la configuración se pueden cambiar usando los elementos de la barra de control. Si desea modificar el método, haga clic en la pestaña **Editar método**.

Con el botón secundario del ratón haga clic en el pocillo y se mostrará el menú contextual:

SM1_	1	SM1_5	SM1_9	SM1_13
	Resu	men		
C	Deta	lles		
	Edita	ir		
5	Edita	r la configuració	ón cinética	
	Copiar configuración cinética			
	Pega	ir configuración	cinética	
S	Gráfi	co: etiquetas mi	últiples	
	Gráfi	co: cinética		
c	Gráfi	co: cinética de e	tiquetas múltipl	es
	Gráfi	co: curvas FLT		
S	Gráfi	co: Spectra		
	Gráfi	co: series de dilu	ución	
C	Ocul	tar o visualizar s	elección	
	Most	trar u ocultar dis	eño	

Al seleccionar **Resumen** se mostrará la ventana siguiente, que ofrece información detallada de la definición y la configuración del pocillo seleccionado:

- Concentraciones	1
E 🚺 Conc. simple. (UA/mL)	
🚺 Conc. simple media. (UA/mL)	
🔚 🚺 Conc. media. (UA/mL)	
Estadísticas	
🖳 🔝 s - Datos de medición	
🖳 🔟 v - Datos de medición	
🚽 🔝 s - Datos de referencia	
🖳 🔽 v - Datos de referencia	
🚽 🔟 s - Datos de diferencia	
🗤 🔔 v - Datos de diferencia	
s - Reducción del blanco	
v - Reducción del blanco	
s - Conc. simple media. (UA/mL)	
🦾 🔟 v - Conc. simple media. (UA/mL)	
arriba	
izquierda derecha	
abajo	
	10

Haga clic en **Siguiente** en la ventana de diseño de placa y se mostrará el cuadro de diálogo **Guardar como**, donde puede escribir un nombre de archivo y observaciones. Para guardar el archivo, haga clic en el botón **Guardar** pequeño, a la izquierda de la ventana; se puede seguir trabajando en el método o en la zona de trabajo. Haga clic en el botón **Guardar**, a la derecha de la parte inferior de la pantalla, para guardar el archivo y cerrar el Asistente. El programa regresará a la lista de Asistentes.



12.1.5 Resumen de la definición de ELISA cuantitativo en magellan

1. Restar valor del blanco

Definiciones en magellan

Haga clic en **Añadir nueva transformación** en la barra de control y se mostrará una ventana para preguntarle si desea definir una **Reducción del blanco**. Haga clic en **S**í y la fórmula de **Reducción del blanco** se asignará automáticamente a todos los pocillos.

2. Definir concentraciones

Definiciones en magellan (Barra de control – Diseño de método / Valores de Conc., Dil. y Ref.)

Identificador seleccionado: ST Unidad: UA/mL

ST1_1	5	(ST1_1Estándar 1 primer grupo experimental)
ST1_2	10	(ST1_2Estándar 2 primer grupo experimental)
ST1_3	20	(ST1_3Estándar 3 primer grupo experimental)
ST1_4	40	(ST1_4Estándar 4 primer grupo experimental)
ST1_5	80	(ST1_5Estándar 5 primer grupo experimental)
ST1_6	160	(ST1_6Estándar 6 primer grupo experimental)

3. Definir curva estándar

Definiciones en magellan (Barra de control – Concentraciones / Curva estándar)

Datos de entrada	reducción del blanco
Tipo de análisis	regresión lineal
Eje X	lineal
eje Y	lineal

4. Definir cortes

Definiciones en magellan (Barra de control – Evaluar datos / Definición de corte)

Datos de entrada: Conc. media (UA/mL)

Límites 22 18

Positivo ≥ 22>intermedio>=18>negativo

N.º de prueba no competitiva

5. Validación QC

Definiciones en magellan (Barra de control – Evaluar datos / Validación QC):

Datos de entrada: Conc. simple (UA/ml)

Condición de validación 1 NC1_1<8 Condición de validación 2 NC1_2<8

NC1_1..... Control negativo primera réplica primer grupo experimental

NC1_2..... Control negativo segunda réplica primer grupo experimental

13. Glosario

Término	Definición
Conc. simple media (???)	Concentración calculada haciendo un promedio de las concentraciones individuales
Base DO	Parámetro cinético: valor base OD/RFU/RLU del cálculo de iniciación
Base DO %	Parámetro cinético: valor base del cálculo de iniciación en %
Portapapeles	El portapapeles es el medio que utilizan los programas de Windows para intercambiar datos entre sí. Se pueden cortar o copiar datos al portapapeles para seleccionarlos en una aplicación de Windows y luego se pueden pegar desde el portapapeles para añadirlos a otra aplicación.
Coef. correlación	El coeficiente de correlación indica la consistencia y la dirección de una relación lineal entre dos variables arbitrarias.
Límite de corte	El límite de corte permite al usuario definir los límites entre dos condiciones (por ejemplo: positivo o intermedio). Estos criterios se utilizan en la evaluación de los resultados.
Resultados de corte	Nombre del rango en el que se encuentran los valores, tal como se haya definido en la definición del corte
Definición del corte	Definición de todos los rangos corte y límites calculados
Factores de dilución	Factores de dilución definidos de las muestras y controles.
	Un factor de dilución de 2 representa una dilución 1:2.
Series de dilución	Una muestra que tiene réplicas con distintas diluciones.
Calidad del ajuste	Uno menos las desviaciones cuadráticas relativas promedio de los puntos base desde la curva
Gráfico	Se pueden mostrar gráficos para mediciones cinéticas, cinéticas de enzimas, mediciones de etiquetas múltiples, series de dilución o curvas estándar.
Oculto	Se muestra si los datos del pocillo están ocultos al imprimir
HUID	Número de identificación de la unidad hardware (Hardware Unit Identification)
IC 50	Dilución/concentración que causa el 50% de la respuesta máxima
No válido	Si un valor es no válido, no admite ningún cálculo
Gráfico: Cinética	Gráfico de mediciones cinéticas
Lámpara tenue	No se obtienen valores medición debido al funcionamiento erróneo del instrumento de absorbencia.
Diseño, Diseño de placa	Define dónde están colocadas en la microplaca las muestras o los controles
Pendiente máx. DO /h	Parámetro cinético: Pendientes máximas de la curva cinética por hora
Pendiente máx. DO /m	Parámetro cinético: Pendientes máximas de la curva cinética por minuto
Pendiente máx. DO /s	Parámetro cinético: Pendientes máximas de la curva cinética por segundo
Máximo DO	Parámetro cinético: Valor máximo de las curvas cinéticas
Pendiente media DO /h	Parámetro cinético: Pendientes medias de la curva cinética por hora



Término	Definición
Pendiente media DO /m	Parámetro cinético: Pendientes medias de la curva cinética por minuto
Pendiente media DO /s	Parámetro cinético: Pendientes medias de la curva cinética por segundo
Conc. media (???)	Concentración calculada a partir de la media de las réplicas de los datos de entrada
Datos de medición	Medición de absorbencia de longitud de onda doble: datos medidos con el filtro de medición
Parámetros de medición	Define el modo de medición, longitud de onda, tamaño de la placa, agitación, etc.
Tipo de medición	El tipo de medición puede ser una medición de punto final, una medición cinética, una medición de etiquetas múltiples o una medición de cinética de pocillo.
Método	Los métodos contienen los parámetros de medición y la definición de la evaluación. Al ejecutar un método se crea una zona de trabajo que contiene los datos medidos y calculados.
Mínimo ???	Parámetro cinético: Valor mínimo de las curvas cinéticas
MultPt	La curva estándar no es monótona y devuelve más de una concentración para los datos de entrada dados.
NoCalc	El cálculo no devuelve ningún valor.
Concentraciones originales	Concentraciones de la curva estándar definida en el método.
Estado de pipeteado	Si se importa una lista de IDs de muestras desde un software de pipeteado, se puede mostrar el estado de pipeteado.
Validación QC	Los criterios de validación QC se definen en el método y estipulan si una medición es válida o no válida. Los criterios pueden, por ejemplo, indicar si los valores de la medición se encuentran demasiado separados entre sí, o si se desvían demasiado de un valor previsto. El programa advierte automáticamente al usuario si no se cumplen los criterios.
Datos sin procesar	Datos medidos por el instrumento.
Datos de referencia	Medición de absorbencia de longitud de onda doble: datos medidos con el filtro de referencia
Estadística de resultados	Resumen del número de valores en los distintos rangos de corte
Lista de IDs de muestras	Los IDs de muestras se asignan a cada pocillo con la idea de que se pueda identificar la sonda asociada. Los IDs por lo general son códigos de barras importados desde listas de IDs de muestras guardadas por un software de pipeteado.
IDs de muestras	Son las identificaciones de las muestras
Conc. simple (???)	Concentración calculada a partir de los datos de entrada de cada réplica
Prueba	En las versiones anteriores, una prueba contenía los ajustes de la evaluación, pero no los parámetros de medición. magellan ya no admite las pruebas, que han sido reemplazadas por métodos más potentes.
Tiempo base DO	Parámetro cinético: tiempo hasta que se alcanza el valor base
Tiempo base DO %	Parámetro cinético: tiempo hasta que se alcanza el % del valor base



Término	Definición
Tiempo base hasta iniciación DO	Parámetro cinético: tiempo entre valor base y valor de iniciación
Tiempo base hasta iniciación DO %	Parámetro cinético: tiempo entre la base y el % del valor de iniciación
Tiempo de pendiente máx. (s)	Parámetro cinético: punto de tiempo de la pendiente máxima
Tiempo de máximo DO	Parámetro cinético: tiempo hasta que se alcanza el máximo
Tiempo de mínimo DO	Parámetro cinético: tiempo hasta que se alcanza el mínimo
Tiempo de iniciación DO	Parámetro cinético: tiempo hasta que se alcanza el valor de iniciación definido
Tiempo de iniciación DO %	Parámetro cinético: tiempo hasta que se alcanza el % de iniciación definido
Puntos de tiempo	Registros de tiempo de las mediciones individuales de una medición cinética
No disponible	Los datos solicitados no están disponibles
No usado	No se ha medido ningún dato, no se ha definido ningún diseño para este pocillo
Indicaciones al usuario	Las indicaciones al usuario son referencias que se configuran en el método. Se muestran antes de la medición y el usuario tiene que aportar texto a las mismas. Ese texto se incorporará más tarde a la copia impresa.
Criterios de validación	Resumen de resultados de condiciones de validación
Posiciones de los pocillos	Nombre del pocillo, por ejemplo: A1, A2,
Zona de trabajo	Todos los datos disponibles dentro del programa se pueden encontrar en la zona de trabajo, por ejemplo, los datos de medición, la definición de impresión y la definición del método. La zona de trabajo se utiliza para cargar métodos y ejecutar mediciones.
!	Precede a los valores eliminados durante el cálculo
#	Precede a los valores de concentración calculados mediante extrapolación y que se encuentran fuera del rango de la curva estándar
()	Los paréntesis rodean los valores que han sido enmascarados
*	Un asterisco marca los valores medidos con la opción "Usar regulación de ganancia", que corrige (=reduce) la ganancia.
??? – Media	Promedios calculados (por ejemplo: Datos sin procesar – Media)
??? – Desviación estándar s - ???	Desviaciones estándar calculadas (por ejemplo: Datos sin procesar – Desviación estándar o s – Datos sin procesar)
??? – Coeficiente de variación v - ???	Coeficientes de variación calculados (por ejemplo: Datos sin procesar – coeficiente de Variación o v – Datos sin procesar)
~	Se añade a valores editados o simulados
<mín< th=""><th>La concentración calculada es inferior al mínimo</th></mín<>	La concentración calculada es inferior al mínimo
>Máx	La concentración calculada supera el máximo
<en blanco=""></en>	Informe impreso: insertar matriz vacía o copia impresa de tabla



Término	Definición
<salto de="" página=""></salto>	Informe impreso: imprimir el siguiente elemento en la página siguiente
<separador></separador>	Informe impreso: imprimir una línea entre dos elementos
x	El símbolo \mathbf{x} se refiere al valor actual dentro de un pocillo
concX	El símbolo concX se refiere a la concentración del estándar en el pocillo actual
dilX	El símbolo dilX se refiere a la dilución de la muestra o el control en el pocillo actual
'???'!	Si hay más de un conjunto de datos de entrada, se utilizarán para los cálculos los datos disponibles
[]	Los índices tienen acceso a los distintos ciclos de una medición cinética, mientras que [0] indica el primer ciclo
*	El asterisco marca los identificadores que se han definido como alias.



Índice

Α

abs(argumento)	181
Acerca de magellan	170
Activar favorito	105
Administración de usuarios	155
Opciones	160
Administración de usuarios (magellan Stand	dard) 163
Administración de usuarios (magellan Track	ker) 155
Cambiar usuario	
Inicio de sesión	164
Administración de usuarios Registro de	
modificaciones	159
Akima	194
algoritmo de Levenberg-Marquardt	198
Añadir HUIDs	173
Añadir nueva transformación	53
Añadir/modificar grupo	158
Añadir/modificar usuario (magellan Standar	d) 164
Añadir/modificar usuario (magellan Tracker)) 156
and – expresión lógica	
Anexar firma	139
Aplicación bloqueada	24
Aprobación	140
Archivado automático	143
Archivo ASTM	81, 82, 83
área bajo la curva cinética	
Asignar alias	40, 45
Asistente	
Elementos estándar	21
Asistente para registrar	170
avg(argumento)	182

В

Bloqueo automático	160
Botón Ayuda	22

С

calcAlways(argumento)	188
Cálculo ICx	49
Cálculos	175
Reducción de datos de espectros	176
Cálculos	175
Cambiar contraseña	24
Caracteres especiales	125
Cinco parámetros – Marquardt	198
Cinética de ciclo arbitraria	109
Cinética de enzima	60
coeficiente de correlación	198
Compatibilidad con lectores	11
Componentes y términos	19
Comprobación de la instalación	13
Concentraciones	127
Conectar un instrumento	25
Configuración de Exportación ASTM	83
Configuración de listas de IDs de muestras .	89
Configurar impresora	120
Confirmación	140
Contraseña	
Opciones	161
Contraseña inicial	24
Control de temperatura1	9, 31, 110
Control del instrumento	31
Convertir a	145
Convertir de	146
Convertir documentos	145

Copiar a Excel	124
Copiar en formato ASCII	124
Correo de prueba	
Corte	66
countDeleted(arg1; arg2)	
Crear o editar IDs de muestras	89
Crear o editar método	33
Criterios de selección de archivos	116
Cuadro de diálogo de bienvenida	22
Cuadro de diálogo Espectros	
Cuatro parámetros	197
Cuatro parámetros – Marquardt	197
Curva estándar	60, 128

D

Datos del instrumento	126
Datos disponibles	
Datos reducidos	126
Datos transformados	126
Definición de delimitador ASTM	81
Definir constantes	
Definir evaluación	
Definir placas de filtro	
Definir un diseño de placas	
Definir un nuevo identificador	
Derechos del usuario	165, 166
Directorios predeterminados	
diseño de método	
Diseño de placas	
1	

Ε

Editar indicaciones al usuario 108 Editar pocillo 93 Ejemplo de aplicación 201 Ejemplo ELISA cuantitativo 201 Elementos estándar 21 eliminate (arg1; arg2; arg3; arg4) 184 eliminatePerc (arg1; arg2; arg3; arg4) 185 esiminateRange (arg1; arg2; arg3; arg4) 185 escala de colores 132 espectro de una medición de exploración 136 Estado de la medición 113 Evaluar datos 66 Evaluar Resultados 115 Evaluar Resultados 138 exp(argumento) 181 Exportar 72 a Excel 73 Exportar a archivo ASCII 120 Exportar a archivo ASTM 120 Exportar a archivo ASTM 120 Exportar a archivo ASTM (LIS) 120 Exportar a LIS 81 expresiones lógicas 180 or 180 ff 180	Editar constantes	108
Editar pocillo 93 Ejemplo de aplicación 201 Ejemplo ELISA cuantitativo 201 Elementos estándar 21 eliminate (arg1; arg2; arg3; arg4) 184 eliminatePerc (arg1; arg2; arg3; arg4) 185 escala de colores 132 espectro de una medición de exploración 136 Estado de la medición 113 Evaluar datos 66 Evaluar Resultados 115 Evaluar Resultados 138 exp(argumento) 181 Exportar a ASCII a ASCII 72 a Excel 73 Exportar a archivo ASTM 120 Exportar a archivo ASTM 120 Exportar a archivo ASTM 120 Exportar a Excel 120 Exportar a LIS 81 expresiones lógicas 180 or 180 F 180	Editar indicaciones al usuario	108
Ejemplo de aplicación 201 Ejemplo ELISA cuantitativo 201 Elementos estándar 21 eliminate (arg1; arg2; arg3; arg4) 184 eliminatePerc (arg1; arg2; arg3; arg4) 185 eliminateRange (arg1; arg2; arg3; arg4) 185 Escala de colores 132 espectro de una medición de exploración 136 Estado de la medición 113 Evaluar datos 66 Evaluar resultados 115 Evaluar Resultados 138 exp(argumento) 181 Exportar 69 a ASCII 72 a Excel 73 Exportar a archivo ASTM 120 Exportar a archivo ASTM 120 Exportar a archivo ASTM 120 Exportar a Excel 120 Exportar a LIS 81 expresiones lógicas 180 or 180 F 180	Editar pocillo	
Ejemplo ELISA cuantitativo 201 Elementos estándar 21 eliminate (arg1; arg2; arg3; arg4) 184 eliminatePerc (arg1; arg2; arg3; arg4) 185 eliminateRange (arg1; arg2; arg3; arg4) 185 escala de colores 132 espectro de una medición de exploración 136 Estado de la medición 113 Evaluar datos 66 Evaluar resultados 115 Evaluar Resultados 138 exp(argumento) 181 Exportar 72 a ASCII 72 a Excel 73 Exportar a archivo ASTM 120 Exportar a archivo ASTM 120 Exportar a archivo ASTM (LIS) 120 Exportar a LIS 81 expresiones lógicas 180 or 180 F 180	Ejemplo de aplicación	201
Elementos estándar. 21 eliminate (arg1; arg2; arg3; arg4) 184 eliminatePerc (arg1; arg2; arg3; arg4) 185 eliminateRange (arg1; arg2; arg3; arg4) 185 Escala de colores 132 espectro de una medición de exploración 136 Estado de la medición 113 Evaluar datos 66 Evaluar resultados 115 Evaluar Resultados 138 exp(argumento) 181 Exportar 72 a ASCII 72 a Excel 73 Exportar a archivo ASTM 120 Exportar a archivo ASTM 120 Exportar a archivo ASTM 120 Exportar a Excel 120 Exportar a archivo ASTM 120 Exportar a archivo ASTM 120 Exportar a LIS 81 Expresiones lógicas 180 or 180 F 180	Ejemplo ELISA cuantitativo	201
eliminate (arg1; arg2; arg3; arg4) 184 eliminatePerc (arg1; arg2; arg3; arg4) 185 eliminateRange (arg1; arg2; arg3; arg4) 185 Escala de colores 132 espectro de una medición de exploración 136 Estado de la medición 113 Evaluar datos 66 Evaluar resultados 115 Evaluar Resultados 138 exp(argumento) 181 Exportaría 69 exportar 72 a ASCII 72 a Excel 73 Exportar a archivo ASTM 120 Exportar a archivo ASTM 120 Exportar a archivo ASTM 120 Exportar a archivo ASTM (LIS) 120 Exportar a LIS 81 expresiones lógicas 180 ad 180 F 180 F Fabricante	Elementos estándar	
eliminatePerc (arg1; arg2; arg3; arg4) 185 eliminateRange (arg1; arg2; arg3; arg4) 185 Escala de colores 132 espectro de una medición de exploración 136 Estado de la medición 113 Evaluar datos 66 Evaluar resultados 115 Evaluar Resultados 138 exp(argumento) 181 Exportación de datos 69 Exportar 72 a ASCII 72 a Excel 73 Exportar a archivo ASTM 120 Exportar a Bxcel 120 Exportar a LIS 81 expresiones lógicas 180 or 180 F 180	eliminate (arg1; arg2; arg3; arg4)	184
eliminateRange (arg1; arg2; arg3; arg4) 185 Escala de colores 132 espectro de una medición de exploración 136 Estado de la medición 113 Evaluar datos 66 Evaluar resultados 115 Evaluar Resultados 138 exp(argumento) 181 Exportación de datos 69 Exportar 72 a ASCII 72 a Excel 73 Exportar a archivo ASCII 120 Exportar a archivo ASTM 120 Exportar a llS 81 Expresiones lógicas 180 or 180 F Fabricante 2	eliminatePerc (arg1; arg2; arg3; arg4)	185
Escala de colores	eliminateRange (arg1; arg2; arg3; arg4)	185
espectro de una medición de exploración 136 Estado de la medición 113 Evaluar datos 66 Evaluar resultados 115 Evaluar Resultados 138 exp(argumento) 181 Exportación de datos 69 Exportar 72 a ASCII 72 a Excel 73 Exportar a archivo ASCII 120 Exportar a archivo ASTM 120 Exportar a bigicas 120 and 180 or 180 F 180 Fabricante 2	Escala de colores	132
Estado de la medición 113 Evaluar datos 66 Evaluar resultados 115 Evaluar Resultados 138 exp(argumento) 181 Exportación de datos 69 Exportar 72 a ASCII 72 a Excel 73 Exportar a archivo ASCII 120 Exportar a archivo ASTM 120 Exportar a llS 81 Expresiones lógicas 180 or 180 F Fabricante	espectro de una medición de exploración	136
Evaluar datos	Estado de la medición	113
Evaluar resultados 115 Evaluar Resultados 138 exp(argumento) 181 Exportación de datos 69 Exportar 69 a ASCII 72 a Excel 73 Exportar a archivo ASCII 120 Exportar a archivo ASTM 120 Exportar a Bit 120 Exportar a LIS 120 Exportar a LIS 81 Expresiones lógicas 180 or 180 ff 180 F Fabricante	Evaluar datos	66
Evaluar Resultados 138 exp(argumento) 181 Exportación de datos 69 Exportar 3 a ASCII 72 a Excel 73 Exportar a archivo ASCII 120 Exportar a archivo ASTM 120 Exportar a archivo ASTM (LIS) 120 Exportar a Excel 120 Exportar a LIS 81 Expresiones lógicas and and 180 or 180 F Fabricante	Evaluar resultados	115
exp(argumento) 181 Exportación de datos 69 Exportar 72 a ASCII 72 a Excel 73 Exportar a archivo ASCII 120 Exportar a archivo ASTM 120 Exportar a archivo ASTM (LIS) 120 Exportar a Excel 120 Exportar a LIS 81 Expresiones lógicas and and 180 F Fabricante	Evaluar Resultados	138
Exportación de datos 69 Exportar 72 a ASCII 72 a Excel 73 Exportar a archivo ASCII 120 Exportar a archivo ASTM 120 Exportar a archivo ASTM (LIS) 120 Exportar a Excel 120 Exportar a LIS 81 Expresiones lógicas 81 and 180 or 180 F Fabricante Fabricante 2	exp(argumento)	181
Exportar a ASCII 72 a Excel 73 Exportar a archivo ASCII 120 Exportar a archivo ASTM 120 Exportar a Excel 120 Exportar a Excel 120 Exportar a LIS 81 Expresiones lógicas and and 180 or 180 F Fabricante	Exportación de datos	69
a ASCII 72 a Excel 73 Exportar a archivo ASCII 120 Exportar a archivo ASTM 120 Exportar a Excel 120 Exportar a Excel 120 Exportar a LIS 81 Expresiones lógicas 81 and 180 or 180 F 180 Fabricante 2	Exportar	
a Excel	a ASCII	72
Exportar a archivo ASCII 120 Exportar a archivo ASTM 120 Exportar a archivo ASTM 120 Exportar a archivo ASTM 120 Exportar a archivo ASTM (LIS) 120 Exportar a Excel 120 Exportar a LIS 81 Expresiones lógicas 180 or 180 F Fabricante Fabricante 2	a Excel	73
Exportar a archivo ASTM	Exportar a archivo ASCII	120
Exportar a archivo ASTM	Exportar a archivo ASTM	120
Exportar a archivo ASTM (LIS)	Exportar a archivo ASTM79, 82, 1	19, 120
Exportar a Excel	Exportar a archivo ASTM (LIS)	120
Exportar a LIS	Exportar a Excel	120
Expresiones lógicas and	Exportar a LIS	
and	Expresiones lógicas	
if() then() else()	and	180
or	if() then() else()	
F Fabricante2	or	
Fabricante2	F	
	Fabricante	2

Aprobación	140
Confirmación	140
Firmar un archivo	139
Formato de números	84
Fórmula	178
Cómo escribir una fórmula	178
Expresiones lógicas	
Funciones	
Funciones básicas	
Funciones de eliminación	
Funciones de espectros	
Funciones estadísticas	
Operadores	
Otras funciones	
Variables	178
Formulario de registro	
frac(argumento)	181
Fuente de impresión	121

G

Glosario	221
Gráfico de cinética	134
Gráfico de cinética de enzima	127
Gráfico de curva estándar	130
Gráfico de etiquetas múltiples	134
Gráfico Levy-Jennings	121
Guardar en archivador	141
Guardar listas de IDs de muestras	104
Guardar los Resultados evaluados	138
Guardar métodos	85

Н

3

I

Icono Otras opciones	141
Opciones	148
Iconos	18
identificadores	45, 46
IDs de muestras	131
if() then() else() – expresión lógica	180
ignore()	187
Importar datos sin procesar	144
Importar lista de IDs de muestras	96
Importar o modificar lista de IDs de muestras.	91
Imprimir	120
In(argumento)	181
Indicaciones al usuario	84
Informe impreso	74
Iniciaciones	58
Iniciar magellan	23
Iniciar medición	105, 108
iniciar sesión	164
Iniciar sesión	24
inicios de sesión	
fallidos	160
Inicios de sesión fallidos	160
Insertar Lista de IDs de muestras	112
Instalación	
Software	12
int(argumento)	181
Interfaz del usuario	17
isInvalid()	187

L

lg(argumento)	181
LIS - Laboratory Information Systems	. 79
Lista de accesos directos	. 22
Lista de asistentes	. 17
lista de IDs de muestras	

importar	. 89	, 96
Log(argumento)	···· [·]	181

Μ

magellan Standard	12
magellan Tracker	12
max(argumento)	183
maxAvg(argumento)	183
mean(argumento)	182
median(argumento)	
medianPlate()	
Medición de etiquetas múltiples	
Mediciones cinéticas	
Mediciones de etiquetas múltiples	
Mediciones de punto final	
Menú contextual de un pocillo en	-
pestaña Evaluar resultados	
Diálogo Detalles	133
Diálogo Gráfico-Cinética	134
Diálogo Gráfico-Espectros	136
Diálogo Gráfico-Etiquetas múltiples	134
Editar diálogo	133
Resumen de un pocilio	40, 132
Menus de la pestana Evaluar resultados	
Metodo de Levenberg-Marquardt	
Método Exportar	
Metodos multiplaca	
min(argumento)	
Min./Máx.	
minAvg(argumento)	183
Movimientos	
portaplacas, placa de filtro	31

Ν

Notas de método	85
-----------------	----

0

Obtener datos sin procesar105,	106
Opciones de Copiar/Exportar	150
Opciones de Correo electrónico	162
or – expresión lógica	180
Organizar favoritos	107

Ρ

Parámetros cinéticos	126
Parametros de medicion	. 34
Pegar desde formato ASCII	124
Pegar en formato ASCII	. 95
Pendientes	. 56
Pestaña Copiar	150
Pestaña Directorios	149
Pestaña Menú contextual de un pocillo de	
Evaluar resultados	132
pestaña Otras opciones	153
Pestaña Otras opciones de Evaluar resultados	132
Pestaña Vista de placa	152
Pestaña Visualizar placa	152
PointwiseCV(argumento)	183
Ponderación	
aiuste de cuatro/cinco parámetros –	
ajuste polinomial/Marguardt	199
Presentación preliminar	120
Procesamiento automatizado de datos	. 79
Procesamiento de archivos	172
Procesamiento de carpetas	20
Procesamiento de datos	69
	. 55
Q	

QC de placa a placa.....69, 121

	•	



R

	•••
Reducción de datos cinéticos	56
Reducción de datos de espectros 50, 1	76
Registrar magellan	27
Registro de modificaciones1	71
Registro de modificaciones del sistema1	73
Reglas Westgard [®] 12	23
Requisitos del sistema	11
Resultados cualitativos1	31
Resumen de administración de usuarios1	62
Resumen de pocillos	95
round(argumento)1	82

S

Seguridad cibernética	
Selección de llenado automático	94
Series de dilución	131
Servidor SMTP	162
símbolos de tipo de archivo	
Software	
Eliminación	15
Instalación	12
Spline cúbica	194

sqr(argumento)	. 182
sort(argumento)	. 182
stddev(argumento)	. 184
Sum(argumento)	. 184

Т

Tipos de análisis de la curva estándar	
Tipos de archivo usados con magellan	
Tipos de mediciones	35
Transformación	
Añadir nueva transformación	53
Cambiar nombre de transformación	53
Datos transformados	53
Transformaciones cinéticas	60, 66
U	
Usar método predefinido	. 105, 107
V	
Validación de operatividad. OQ	13. 16
Validación OC	68 132
	00, 102

Validación QC	68, 132
Valores Conc., Dil. y Ref	
Ventana de diseño de placas	40, 125
Vista de conjunto de la zona de trabajo	
Volver a calcular con otro método	

Marcas registradas

Los siguientes nombres de productos y todas las marcas comerciales registradas o no que se mencionan en este documento se utilizan únicamente a efectos de identificación y son propiedad exclusiva de sus respectivos titulares:

- [Nombre de producto][™] (infinite®) Tecan® y el logotipo de Tecan son marcas registradas de Tecan Group Ltd., Männedorf, Switzerland
- DNA Expert[™] es una marca comercial registrada de Techcomp Ltd., Hong Kong, China
- Agilent® es una marca comercial registrada de Agilent Technologies, Inc., Santa Clara, CA, USA
- AIR LIQUIDE[™] es una marca comercial registrada de AIR LIQUIDE, S.A., Paris, France
- Aseptisol® es una marca comercial registrada de BODE Chemie GmbH & Co. KG, Hamburg, Germany
- Bacillol® es una marca comercial registrada de BODE Chemie GmbH & Co. KG, Hamburg, Germany
- Costar®, Corning® y NBS[™] son marcas registradas de Corning Incorporated, Corning, NY, USA
- Greiner®, µClear®, Lumitrac[™] y Fluotrac[™] son marcas registradas de Greiner Labortechnik GmbH, Frickenhausen, Germany
- HTRF® es una marca comercial registrada de Cisbio International, France
- Invitrogen[™] es una marca comercial registrada de Invitrogen Corporation, Carlsbad, CA, USA
- Lysetol® y Gigasept® (formerly Lysetol) son marcas registradas de Schülke & Mayr GmbH, Norderstedt, Germany
- Microcide® es una marca comercial registrada de Global Biotechnologies Inc., Portland, Maine, USA
- Microman® es una marca comercial registrada de Gilson, Inc., Middleton, WI, USA
- Pentium® es una marca comercial registrada de Intel Corporation, Santa Clara, CA, USA
- Invitrogen® y PanVera® son marcas registradas de Invitrogen Corporation Carlsbad, CA, USA
- Windows®, MS DOS®, Visual Basic® y Excel® son marcas registradas de Microsoft Corporation, Redmond, WA, USA
- Hamamatsu® es una marca comercial registrada de HAMAMATSU Photonics K.K. [IR], Hamamatsu City, Japan
- NUNC[™] y Matrix son marcas registradas de Thermo Fisher Scientific, Waltham, MA, USA
- Polyfiltronics® es una marca comercial registrada de Whatman International Ltd.
- Dynex® es una marca comercial registrada de Magellan Biosciences, Chelmsford, MA, USA
- Labsystem® es una marca comercial registrada de Labsystem kft., Budapest, Hungary
- BRET²®, DeepBlueC® y PerkinElmer® son marcas registradas de PerkinElmer, Inc., Waltham, Massachusetts, USA
- Chroma-Glo[™] es una marca comercial registrada de Promega Corporation, Madison, WI 53711 USA
- MycoAlert® es una marca comercial registrada de Cambrex Corporation, East Rutherford, NJ, USA

Atención al Cliente de Tecan

Si necesita soporte técnico o tiene consultas relativas a su producto de Tecan, póngase en contacto con su servicio local de Atención al Cliente de Tecan. Consulte la información de contacto en <u>http://www.tecan.com/</u>.

Antes de ponerse en contacto con Tecan para solicitar soporte técnico, prepare la siguiente información para obtener un soporte óptimo (consulte la placa de características):

- Modelo de su producto
- Número de serie (SN) de su producto
- Software y versión del software (cuando corresponda)
- Descripción del problema y persona de contacto
- Fecha y hora de aparición del problema
- Medidas que ya ha tomado para corregir el problema
- Su información de contacto (teléfono, fax, correo electrónico, etc.)

End User Software License Agreement

IMPORTANT-READ CAREFULLY: This End-User License Agreement ("EULA") is a legal agreement between you (either an individual or a legal entity) and Tecan Austria GmbH ("Tecan") for the proprietary software product identified above, which includes computer software ("SOFTWARE PRODUCT") and may include associated media, printed materials, and "online" or electronic documentation ("DOCUMENTATION"). This SOFTWARE PRODUCT is made available to you only on the terms and conditions of this EULA. By installing, copying, or otherwise using the SOFTWARE PRODUCT (which ever occurs first), you agree to be bound by the terms of this EULA. If you do not agree with the terms of this EULA, you are not authorized to install and/or use the SOFTWARE PRODUCT.

SOFTWARE PRODUCT LICENSE

The SOFTWARE PRODUCT and the DOCUMENTATION are protected by copyright laws and international copyright treaties, as well as other intellectual property laws and treaties. The SOFTWARE PRODUCT is licensed, not sold.

1. GRANT OF LICENSE

This EULA grants you the non-exclusive and non-transferable right to use the SOFTWARE PRODUCT in accordance with the instructions and procedures in the DOCUMENTATION for your own internal purposes only. You may install and use the number of copies of the SOFTWARE PRODUCT on the number of computers or workstations connected to the number of devices as specified in the documentation to this EULA in accordance with the software installation procedure described in the documentation.

You may either make one copy of the SOFTWARE PRODUCT solely for archival purposes in support of your use of the SOFTWARE PRODUCT on a single computer or transfer the SOFTWARE PRODUCT to a single hard disk provided that you keep the original only for backup or archival purposes. You may not reproduce or distribute the SOFTWARE PRODUCT in any other way, including rental or leasing.

The SOFTWARE PRODUCT shall be used exclusively with the devices specified in the documentation and must not be used on any other devices.

2. LICENSE FEE

The license fee for the right to use the SOFTWARE PRODUCT is set forth in the documentation.

3. DESCRIPTION OF OTHER RIGHTS AND LIMITATIONS

You may not reverse engineer, decompile, or disassemble the SOFTWARE PRODUCT, except and only to the extent that such activity is expressly permitted by applicable law notwithstanding this limitation.

The SOFTWARE PRODUCT is licensed as a single product. Its component parts may not be separated for use on more than one computer or workstation.

Tecan may provide you with support services related to the SOFTWARE PRODUCT ("SUPPORT SERVICES"). Use of SUPPORT SERVICES is governed by Tecan's policies and programs described in the user manual, in "online" documentation, and/or in other materials provided by Tecan. Any supplemental software code provided to you as part of the SUPPORT SERVICES shall be considered part of the SOFTWARE PRODUCT and subject to the terms and conditions of this EULA. With respect to technical information you provide to Tecan as part of the SUPPORT SERVICES, Tecan may use such information for its business purposes, including for product support and development. Tecan will not utilize such technical information in a form that personally identifies you.

You may not transfer any of your rights under this EULA without the prior written consent of Tecan to a third party.

Without prejudice to any other rights, Tecan may terminate this EULA if you fail to comply with the terms and conditions of this EULA. In such event, you may not use the SOFTWARE PRODUCT any longer and you must destroy all copies of the SOFTWARE PRODUCT and the DOCUMENTATION and all of its component parts within a period of fourteen days.

4. COPYRIGHT

All title and copyrights in and to the SOFTWARE PRODUCT (including but not limited to any charts, images, photographs, animations, video, audio, music, text, and "applets" incorporated into the SOFTWARE PRODUCT), the DOCUMENTATION and any copies of the SOFTWARE PRODUCT and the DOCUMENTATION are owned by Tecan Austria GmbH or its suppliers and licensed to Tecan. The SOFTWARE PRODUCT and the DOCUMENTATION are protected by copyright laws and international treaty provisions. Therefore, you must treat the SOFTWARE PRODUCT and the DOCUMENTATION are the DOCUMENTATION like any other copyright protected material.

5. TERMS AND CONDITIONS FOR USE

Use of the SOFTWARE PRODUCT shall be subject to compliance with the following terms and conditions relating thereto:

- The SOFTWARE PRODUCT shall be used exclusively on devices specified in the documentation;

- only the latest version of the SOFTWARE PRODUCT shall be used;

- the SOFTWARE PRODUCT shall not be modified nor caused to be modified.

It shall be your responsibility to personally effect any requisite declarations to the authorities with a view to using SOFTWARE PRODUCTS.



6. PRECAUTIONS OF USE

Before being commercialized, the SOFTWARE PRODUCT underwent a series of tests to measure the reliability of results obtained.

However, taking into account the high number of possible applications for which the SOFTWARE PRODUCT can be used, it has not been possible to carry out said tests in a real work situation.

The contracting parties agree that it is not possible in practice to produce data processing programs which will be 100% suitable for all applications.

Therefore, you are strongly advised to verify and to validate results obtained before using the SOFTWARE PRODUCT in a real work situation every time it is used in a new application.

Should there be any doubt as to the results, or in the case of erroneous results, you are requested to immediately contact Tecan.

For the perfect use of SOFTWARE PRODUCTS you are reminded that:

- You must have the requisite knowledge;

- you must, in a regular fashion, write-protect and backup, in particular, to protect the data files from incidents such as deletion, overwriting, virus infection, etc.

7. LIMITED WARRANTY

The SOFTWARE PRODUCT is under warranty for a period of 90 days as of the date of delivery thereof.

During said period, Tecan shall use reasonable efforts to deliver, as quickly as possible and by any means it chooses an updated version or a copy of the said version of the program free of any significant defect appeared.

The provisions mentioned above determine the extent of the warranty granted to you.

Note on Java Support. The SOFTWARE PRODUCT may contain support for programs written in Java. Java technology is not fault tolerant and is not designed, manufactured, or intended for use or resale as online control equipment in hazardous environments requiring fail-safe performance, such as in the operation of nuclear facilities, aircraft navigation or communication systems, air traffic control, direct life support machines, or weapons systems, in which the failure of Java technology could lead directly to death, personal injury, or severe physical or environmental damage. Sun Microsystems, Inc. requires the addition of this disclaimer.

8. RESTRICTED LIABILITY

Tecan shall be subject to an undertaking to exercise due-care.

Said undertaking shall concern the supply of the software conforming to the functionalities described in the documentation thereof, no guarantee whatsoever being given either with regard to results obtained or the fitness of the SOFTWARE PRODUCT for your purposes.

You shall be solely liable for the use of the SOFTWARE PRODUCT and results obtained. In particular, you shall be responsible for verifying the contents and the consistency thereof, as the SOFTWARE PRODUCT shall be used under your sole management, control and responsibility.

TECAN SHALL NOT BE HELD LIABLE FOR ANY INDIRECT OR CONSEQUENTIAL DAMAGES OR ANY LOSS OF PROFIT, SHORTFALL IN EARNINGS OR LOSS OF DATA INCURRED BY YOU EVEN IF TECAN HAS BEEN INFORMED THAT THE SAID LOSS OR PREJUDICE COULD OCCUR.

The SOFTWARE PRODUCT has undergone an anti-virus test. However, Tecan shall not be liable should a virus be present that was undetectable by the anti-virus utility used to run the test mentioned above.

SHOULD FOR ANY REASON WHATSOEVER, TECAN BE HELD LIABLE BY THE COURTS, THE AMOUNT OF ANY DAMAGES EFFECTIVELY OWED AND ORDERED TO BE PAID SHALL, WITH THE EXCEPTION OF INTENTIONAL MISCONDUCT OR GROSS NEGLIGENCE, NOT EXCEED THE TOTAL AMOUNT OF THE SUMS RECEIVED BY TECAN BY VIRTUE OF THE PRESENT EULA.

9. CONFIDENTIALITY

You hereby acknowledge and agree that you will not disclose the SOFTWARE PRODUCT and the DOCUMENTATION to any of your employees (except to those responsible for the application) or to any third party without the prior written consent of Tecan and that this duty of confidentiality survives the term of this EULA. You will safeguard the SOFTWARE PRODUCT and the DOCUMENTATION with the highest degree of care.

10. TERMINATION

If you fail to comply with any of your obligations hereunder, Tecan shall have the right, at any time, to terminate the EULA and take immediate possession of the SOFTWARE PRODUCT and the DOCUMENTATION and of all copies wherever located without demand or notice.

TECAN AUSTRIA GMBH, Untersbergstrasse 1a, A-5082 Grödig / Salzburg, Austria T +43 62 46 89 330, F +43 62 46 72 770, office.austria@tecan.com, www.tecan.com



Declaration of Conformity

We, TECAN Austria GmbH herewith declare under our sole responsibility that the product identified as:

Product Type: Software

Model Designation: SW MAGELLAN V7.5

Article Numbers: 30143623, 30143624, 30143625, 30143626, 30143627, 30143628

Address: Tecan Austria GmbH Untersbergstr. 1A

A-5082 Grödig, Austria SRN: AT-MF-000020241

is in conformity with the provisions of the following EC Directive(s)/Regulation(s) when installed in accordance with the installation instructions contained in the product documentation:

Regulation IVD-R

on in vitro diagnostic devices

For products placed on the Swiss market the CH authorized representative is:

Tecan Schweiz AG Seestrasse 103, 8708 Männedorf, Switzerland



is in conformity with the relevant U.K. legislation for UKCA-marking when installed in accordance with the installation instructions contained in the product documentation:

Medical Devices Regulations 2002

Classification: Other device (all devices except Annex II and self-testing devices) Conformity assessment procedure: Annex III

For products placed on the United Kingdom market the **UK responsible person** is: Tecan UK Limited Theale Court, 11-13 High Street, Theale Reading, Berkshire, RG7 5AH United Kingdom

The current applicable versions of the directives and regulations as well as the list of applied standards which were taken in consideration can be found in separate CE & UK declarations of conformity.

These Instructions for Use and the included Declaration of Conformity are valid for all Magellan SW Packages with the article numbers listed above.